
Orders Report

Here is a simple example that shows how to create a table of contents for your report. There are various ways to build tables of contents. Just let free your imagination and be creative, because JasperReports is very flexible and won't let you down. Most of the time you will end up using scriptlets and subreports, like we have in this sample.

Note, that normally, the table of contents is generated as the report itself is being built. In this sample, it is generated using a subreport placed on the master report summary section and using data prepared by the master report scriptlet object. After the report is generated, the table of contents appears at the end of the master report, as you would expect after checking the report design.

But the parent application performs a trick or a hack, or whatever you want to call it to remove the table of contents from the end of the document and insert it at the beginning.

This operation has apparently only one inconvenience. The final document does not reflect the modification in the order of pages. You can see that in our sample, the table of contents does not display a page number on the footer and that the page #1 comes after the table of contents.

Table of Contents

| | | |
|-----------------|-----------|-----------|
| Letter A | | 1 |
| Argentina | | 1 |
| Austria | | 1 |
| Letter B | | 2 |
| Belgium | | 2 |
| Brazil | | 3 |
| Letter C | | 6 |
| Canada | | 6 |
| Letter D | | 7 |
| Denmark | | 7 |
| Letter F | | 7 |
| Finland | | 7 |
| France | | 8 |
| Letter G | | 10 |
| Germany | | 10 |
| Letter I | | 14 |
| Ireland | | 14 |
| Italy | | 15 |
| Letter M | | 16 |
| Mexico | | 16 |
| Letter N | | 17 |

Table of Contents

| | | |
|-----------------|-----------|-----------|
| Norway | | 17 |
| Letter P | <hr/> | 17 |
| Poland | | 17 |
| Portugal | | 17 |
| Letter S | <hr/> | 18 |
| Spain | | 18 |
| Sweden | | 18 |
| Switzerland | | 20 |
| Letter U | <hr/> | 20 |
| UK | | 20 |
| USA | | 22 |
| Letter V | <hr/> | 26 |
| Venezuela | | 26 |

| ID | Name | City |
|----|------|------|
|----|------|------|

Countries Starting With Letter : A

1. Argentina

| | | |
|-------|----------------------------|--------------|
| 11054 | Cactus Comidas para llevar | Buenos Aires |
| 11019 | Rancho grande | Buenos Aires |
| 10986 | Océano Atlántico Ltda. | Buenos Aires |
| 10958 | Océano Atlántico Ltda. | Buenos Aires |
| 10937 | Cactus Comidas para llevar | Buenos Aires |
| 10916 | Rancho grande | Buenos Aires |
| 10898 | Océano Atlántico Ltda. | Buenos Aires |
| 10881 | Cactus Comidas para llevar | Buenos Aires |
| 10828 | Rancho grande | Buenos Aires |
| 10819 | Cactus Comidas para llevar | Buenos Aires |
| 10782 | Cactus Comidas para llevar | Buenos Aires |
| 10716 | Rancho grande | Buenos Aires |
| 10531 | Océano Atlántico Ltda. | Buenos Aires |
| 10521 | Cactus Comidas para llevar | Buenos Aires |
| 10448 | Rancho grande | Buenos Aires |
| 10409 | Océano Atlántico Ltda. | Buenos Aires |

Count : 16

2. Austria

| | | |
|-------|------------------|----------|
| 11072 | Ernst Handel | Graz |
| 11053 | Piccolo und mehr | Salzburg |
| 11017 | Ernst Handel | Graz |
| 11008 | Ernst Handel | Graz |
| 10990 | Ernst Handel | Graz |
| 10979 | Ernst Handel | Graz |
| 10968 | Ernst Handel | Graz |
| 10895 | Ernst Handel | Graz |
| 10854 | Ernst Handel | Graz |
| 10844 | Piccolo und mehr | Salzburg |
| 10836 | Ernst Handel | Graz |
| 10795 | Ernst Handel | Graz |
| 10776 | Ernst Handel | Graz |

| ID | Name | City |
|-------|------------------|----------|
| 10773 | Ernst Handel | Graz |
| 10771 | Ernst Handel | Graz |
| 10764 | Ernst Handel | Graz |
| 10747 | Piccolo und mehr | Salzburg |
| 10698 | Ernst Handel | Graz |
| 10686 | Piccolo und mehr | Salzburg |
| 10667 | Ernst Handel | Graz |
| 10633 | Ernst Handel | Graz |
| 10597 | Piccolo und mehr | Salzburg |
| 10595 | Ernst Handel | Graz |
| 10571 | Ernst Handel | Graz |
| 10530 | Piccolo und mehr | Salzburg |
| 10514 | Ernst Handel | Graz |
| 10489 | Piccolo und mehr | Salzburg |
| 10442 | Ernst Handel | Graz |
| 10430 | Ernst Handel | Graz |
| 10427 | Piccolo und mehr | Salzburg |
| 10402 | Ernst Handel | Graz |
| 10403 | Ernst Handel | Graz |
| 10392 | Piccolo und mehr | Salzburg |
| 10390 | Ernst Handel | Graz |
| 10382 | Ernst Handel | Graz |
| 10368 | Ernst Handel | Graz |
| 10353 | Piccolo und mehr | Salzburg |
| 10351 | Ernst Handel | Graz |
| 10263 | Ernst Handel | Graz |
| 10258 | Ernst Handel | Graz |

Count : 40

Countries Starting With Letter : B

3. Belgium

| | | |
|-------|------------------|-----------|
| 11038 | Suprêmes délices | Charleroi |
| 11035 | Suprêmes délices | Charleroi |
| 11004 | Maison Dewey | Bruxelles |

| ID | Name | City |
|-------|------------------|-----------|
| 10978 | Maison Dewey | Bruxelles |
| 10930 | Suprêmes délices | Charleroi |
| 10896 | Maison Dewey | Bruxelles |
| 10892 | Maison Dewey | Bruxelles |
| 10885 | Suprêmes délices | Charleroi |
| 10846 | Suprêmes délices | Charleroi |
| 10841 | Suprêmes délices | Charleroi |
| 10767 | Suprêmes délices | Charleroi |
| 10760 | Maison Dewey | Bruxelles |
| 10649 | Maison Dewey | Bruxelles |
| 10529 | Maison Dewey | Bruxelles |
| 10475 | Suprêmes délices | Charleroi |
| 10463 | Suprêmes délices | Charleroi |
| 10458 | Suprêmes délices | Charleroi |
| 10302 | Suprêmes délices | Charleroi |
| 10252 | Suprêmes délices | Charleroi |

Count : 19

4. Brazil

| | | |
|-------|------------------------|----------------|
| 11068 | Queen Cozinha | Sao Paulo |
| 11059 | Ricardo Adocicados | Rio de Janeiro |
| 11052 | Hanari Carnes | Rio de Janeiro |
| 11049 | Gourmet Lanchonetes | Campinas |
| 11042 | Comércio Mineiro | Sao Paulo |
| 11022 | Hanari Carnes | Rio de Janeiro |
| 10989 | Que Delícia | Rio de Janeiro |
| 10981 | Hanari Carnes | Rio de Janeiro |
| 10969 | Comércio Mineiro | Sao Paulo |
| 10961 | Queen Cozinha | Sao Paulo |
| 10959 | Gourmet Lanchonetes | Campinas |
| 10935 | Wellington Importadora | Resende |
| 10925 | Hanari Carnes | Rio de Janeiro |
| 10922 | Hanari Carnes | Rio de Janeiro |
| 10914 | Queen Cozinha | Sao Paulo |
| 10913 | Queen Cozinha | Sao Paulo |

| ID | Name | City |
|-------|------------------------|----------------|
| 10905 | Wellington Importadora | Resende |
| 10903 | Hanari Carnes | Rio de Janeiro |
| 10900 | Wellington Importadora | Resende |
| 10886 | Hanari Carnes | Rio de Janeiro |
| 10877 | Ricardo Adocicados | Rio de Janeiro |
| 10868 | Queen Cozinha | Sao Paulo |
| 10851 | Ricardo Adocicados | Rio de Janeiro |
| 10839 | Tradição Hipermercados | Sao Paulo |
| 10834 | Tradição Hipermercados | Sao Paulo |
| 10830 | Tradição Hipermercados | Sao Paulo |
| 10813 | Ricardo Adocicados | Rio de Janeiro |
| 10809 | Wellington Importadora | Resende |
| 10803 | Wellington Importadora | Resende |
| 10794 | Que Delícia | Rio de Janeiro |
| 10790 | Gourmet Lanchonetes | Campinas |
| 10786 | Queen Cozinha | Sao Paulo |
| 10783 | Hanari Carnes | Rio de Janeiro |
| 10777 | Gourmet Lanchonetes | Campinas |
| 10770 | Hanari Carnes | Rio de Janeiro |
| 10734 | Gourmet Lanchonetes | Campinas |
| 10728 | Queen Cozinha | Sao Paulo |
| 10725 | Familia Arquibaldo | Sao Paulo |
| 10720 | Que Delícia | Rio de Janeiro |
| 10709 | Gourmet Lanchonetes | Campinas |
| 10704 | Queen Cozinha | Sao Paulo |
| 10690 | Hanari Carnes | Rio de Janeiro |
| 10685 | Gourmet Lanchonetes | Campinas |
| 10659 | Queen Cozinha | Sao Paulo |
| 10652 | Gourmet Lanchonetes | Campinas |
| 10650 | Familia Arquibaldo | Sao Paulo |
| 10648 | Ricardo Adocicados | Rio de Janeiro |
| 10647 | Que Delícia | Rio de Janeiro |
| 10644 | Wellington Importadora | Resende |
| 10645 | Hanari Carnes | Rio de Janeiro |

| ID | Name | City |
|-------|------------------------|----------------|
| 10637 | Queen Cozinha | Sao Paulo |
| 10622 | Ricardo Adocicados | Rio de Janeiro |
| 10606 | Tradição Hipermercados | Sao Paulo |
| 10587 | Que Delícia | Rio de Janeiro |
| 10585 | Wellington Importadora | Resende |
| 10581 | Familia Arquibaldo | Sao Paulo |
| 10563 | Ricardo Adocicados | Rio de Janeiro |
| 10541 | Hanari Carnes | Rio de Janeiro |
| 10512 | Familia Arquibaldo | Sao Paulo |
| 10496 | Tradição Hipermercados | Sao Paulo |
| 10494 | Comércio Mineiro | Sao Paulo |
| 10487 | Queen Cozinha | Sao Paulo |
| 10481 | Ricardo Adocicados | Rio de Janeiro |
| 10466 | Comércio Mineiro | Sao Paulo |
| 10447 | Ricardo Adocicados | Rio de Janeiro |
| 10423 | Gourmet Lanchonetes | Campinas |
| 10420 | Wellington Importadora | Resende |
| 10421 | Que Delícia | Rio de Janeiro |
| 10414 | Familia Arquibaldo | Sao Paulo |
| 10406 | Queen Cozinha | Sao Paulo |
| 10386 | Familia Arquibaldo | Sao Paulo |
| 10379 | Que Delícia | Rio de Janeiro |
| 10372 | Queen Cozinha | Sao Paulo |
| 10347 | Familia Arquibaldo | Sao Paulo |
| 10299 | Ricardo Adocicados | Rio de Janeiro |
| 10292 | Tradição Hipermercados | Sao Paulo |
| 10290 | Comércio Mineiro | Sao Paulo |
| 10291 | Que Delícia | Rio de Janeiro |
| 10287 | Ricardo Adocicados | Rio de Janeiro |
| 10261 | Que Delícia | Rio de Janeiro |
| 10256 | Wellington Importadora | Resende |
| 10253 | Hanari Carnes | Rio de Janeiro |
| 10250 | Hanari Carnes | Rio de Janeiro |

Count : 83

| ID | Name | City |
|---|-------------------------------|-----------|
| <i>Countries Starting With Letter : C</i> | | |
| 5. Canada | | |
| 11048 | Bottom-Dollar Markets | Tsawassen |
| 11045 | Bottom-Dollar Markets | Tsawassen |
| 11027 | Bottom-Dollar Markets | Tsawassen |
| 10982 | Bottom-Dollar Markets | Tsawassen |
| 10975 | Bottom-Dollar Markets | Tsawassen |
| 10949 | Bottom-Dollar Markets | Tsawassen |
| 10944 | Bottom-Dollar Markets | Tsawassen |
| 10918 | Bottom-Dollar Markets | Tsawassen |
| 10810 | Laughing Bacchus Wine Cellars | Vancouver |
| 10742 | Bottom-Dollar Markets | Tsawassen |
| 10724 | Mère Paillarde | Montréal |
| 10620 | Laughing Bacchus Wine Cellars | Vancouver |
| 10618 | Mère Paillarde | Montréal |
| 10619 | Mère Paillarde | Montréal |
| 10605 | Mère Paillarde | Montréal |
| 10590 | Mère Paillarde | Montréal |
| 10570 | Mère Paillarde | Montréal |
| 10565 | Mère Paillarde | Montréal |
| 10505 | Mère Paillarde | Montréal |
| 10495 | Laughing Bacchus Wine Cellars | Vancouver |
| 10492 | Bottom-Dollar Markets | Tsawassen |
| 10439 | Mère Paillarde | Montréal |
| 10431 | Bottom-Dollar Markets | Tsawassen |
| 10424 | Mère Paillarde | Montréal |
| 10410 | Bottom-Dollar Markets | Tsawassen |
| 10411 | Bottom-Dollar Markets | Tsawassen |
| 10389 | Bottom-Dollar Markets | Tsawassen |
| 10376 | Mère Paillarde | Montréal |
| 10339 | Mère Paillarde | Montréal |
| 10332 | Mère Paillarde | Montréal |

Count : 30

| ID | Name | City |
|----|------|------|
|----|------|------|

Countries Starting With Letter : D

6. Denmark

| | | |
|-------|---------------|-----------|
| 11074 | Simons bistro | Kobenhavn |
| 10994 | Vaffeljernet | Århus |
| 10946 | Vaffeljernet | Århus |
| 10921 | Vaffeljernet | Århus |
| 10802 | Simons bistro | Kobenhavn |
| 10769 | Vaffeljernet | Århus |
| 10744 | Vaffeljernet | Århus |
| 10688 | Vaffeljernet | Århus |
| 10669 | Simons bistro | Kobenhavn |
| 10642 | Simons bistro | Kobenhavn |
| 10602 | Vaffeljernet | Århus |
| 10591 | Vaffeljernet | Århus |
| 10556 | Simons bistro | Kobenhavn |
| 10465 | Vaffeljernet | Århus |
| 10417 | Simons bistro | Kobenhavn |
| 10399 | Vaffeljernet | Århus |
| 10367 | Vaffeljernet | Århus |
| 10341 | Simons bistro | Kobenhavn |

Count : 18

Countries Starting With Letter : F

7. Finland

| | | |
|-------|----------------|----------|
| 11025 | Wartian Herkku | Oulu |
| 11005 | Wilman Kala | Helsinki |
| 10910 | Wilman Kala | Helsinki |
| 10879 | Wilman Kala | Helsinki |
| 10873 | Wilman Kala | Helsinki |
| 10781 | Wartian Herkku | Oulu |
| 10750 | Wartian Herkku | Oulu |
| 10695 | Wilman Kala | Helsinki |
| 10673 | Wilman Kala | Helsinki |
| 10636 | Wartian Herkku | Oulu |

| ID | Name | City |
|-------|----------------|----------|
| 10615 | Wilman Kala | Helsinki |
| 10583 | Wartian Herkku | Oulu |
| 10553 | Wartian Herkku | Oulu |
| 10526 | Wartian Herkku | Oulu |
| 10455 | Wartian Herkku | Oulu |
| 10437 | Wartian Herkku | Oulu |
| 10416 | Wartian Herkku | Oulu |
| 10412 | Wartian Herkku | Oulu |
| 10333 | Wartian Herkku | Oulu |
| 10320 | Wartian Herkku | Oulu |
| 10270 | Wartian Herkku | Oulu |
| 10266 | Wartian Herkku | Oulu |

Count : 22

8. France

| | | |
|-------|----------------------|------------|
| 11076 | Bon app' | Marseille |
| 11051 | La maison d'Asie | Toulouse |
| 11043 | Spécialités du monde | Paris |
| 10972 | La corne d'abondance | Versailles |
| 10973 | La corne d'abondance | Versailles |
| 10971 | France restauration | Nantes |
| 10964 | Spécialités du monde | Paris |
| 10940 | Bon app' | Marseille |
| 10932 | Bon app' | Marseille |
| 10927 | La corne d'abondance | Versailles |
| 10923 | La maison d'Asie | Toulouse |
| 10907 | Spécialités du monde | Paris |
| 10890 | Du monde entier | Nantes |
| 10876 | Bon app' | Marseille |
| 10871 | Bon app' | Marseille |
| 10860 | France restauration | Nantes |
| 10858 | La corne d'abondance | Versailles |
| 10850 | Victuailles en stock | Lyon |
| 10843 | Victuailles en stock | Lyon |
| 10832 | La maison d'Asie | Toulouse |

| ID | Name | City |
|-------|---------------------------|------------|
| 10826 | Blondel père et fils | Strasbourg |
| 10827 | Bon app' | Marseille |
| 10814 | Victuailles en stock | Lyon |
| 10806 | Victuailles en stock | Lyon |
| 10789 | Folies gourmandes | Lille |
| 10787 | La maison d'Asie | Toulouse |
| 10763 | Folies gourmandes | Lille |
| 10755 | Bon app' | Marseille |
| 10738 | Spécialités du monde | Paris |
| 10739 | Vins et alcools Chevalier | Reims |
| 10737 | Vins et alcools Chevalier | Reims |
| 10732 | Bon app' | Marseille |
| 10730 | Bon app' | Marseille |
| 10715 | Bon app' | Marseille |
| 10683 | Du monde entier | Nantes |
| 10679 | Blondel père et fils | Strasbourg |
| 10671 | France restauration | Nantes |
| 10663 | Bon app' | Marseille |
| 10634 | Folies gourmandes | Lille |
| 10631 | La maison d'Asie | Toulouse |
| 10628 | Blondel père et fils | Strasbourg |
| 10610 | La maison d'Asie | Toulouse |
| 10609 | Du monde entier | Nantes |
| 10584 | Blondel père et fils | Strasbourg |
| 10566 | Blondel père et fils | Strasbourg |
| 10559 | Blondel père et fils | Strasbourg |
| 10546 | Victuailles en stock | Lyon |
| 10525 | Bon app' | Marseille |
| 10511 | Bon app' | Marseille |
| 10500 | La maison d'Asie | Toulouse |
| 10493 | La maison d'Asie | Toulouse |
| 10480 | Folies gourmandes | Lille |
| 10478 | Victuailles en stock | Lyon |
| 10470 | Bon app' | Marseille |

| ID | Name | City |
|-------|---------------------------|------------|
| 10459 | Victuailles en stock | Lyon |
| 10454 | La maison d'Asie | Toulouse |
| 10450 | Victuailles en stock | Lyon |
| 10449 | Blondel père et fils | Strasbourg |
| 10436 | Blondel père et fils | Strasbourg |
| 10425 | La maison d'Asie | Toulouse |
| 10413 | La maison d'Asie | Toulouse |
| 10408 | Folies gourmandes | Lille |
| 10371 | La maison d'Asie | Toulouse |
| 10362 | Bon app' | Marseille |
| 10360 | Blondel père et fils | Strasbourg |
| 10358 | La maison d'Asie | Toulouse |
| 10350 | La maison d'Asie | Toulouse |
| 10340 | Bon app' | Marseille |
| 10334 | Victuailles en stock | Lyon |
| 10331 | Bon app' | Marseille |
| 10311 | Du monde entier | Nantes |
| 10297 | Blondel père et fils | Strasbourg |
| 10295 | Vins et alcools Chevalier | Reims |
| 10274 | Vins et alcools Chevalier | Reims |
| 10265 | Blondel père et fils | Strasbourg |
| 10251 | Victuailles en stock | Lyon |
| 10248 | Vins et alcools Chevalier | Reims |

Count : 77

Countries Starting With Letter : G

9. Germany

| | | |
|-------|--------------------------|----------------|
| 11070 | Lehmanns Marktstand | Frankfurt a.M. |
| 11067 | Drachenblut Delikatessen | Aachen |
| 11058 | Blauer See Delikatessen | Mannheim |
| 11046 | Die Wandernde Kuh | Stuttgart |
| 11036 | Drachenblut Delikatessen | Aachen |
| 11028 | Königlich Essen | Brandenburg |
| 11020 | Ottilies Käseladen | Köln |

| ID | Name | City |
|-------|--------------------------|----------------|
| 11021 | QUICK-Stop | Cunewalde |
| 11012 | Frankenversand | München |
| 11011 | Alfred's Futterkiste | Berlin |
| 10999 | Ottilies Käseladen | Köln |
| 10996 | QUICK-Stop | Cunewalde |
| 10991 | QUICK-Stop | Cunewalde |
| 10967 | Toms Spezialitäten | Münster |
| 10962 | QUICK-Stop | Cunewalde |
| 10956 | Blauer See Delikatessen | Mannheim |
| 10952 | Alfred's Futterkiste | Berlin |
| 10945 | Morgenstern Gesundkost | Leipzig |
| 10938 | QUICK-Stop | Cunewalde |
| 10934 | Lehmanns Marktstand | Frankfurt a.M. |
| 10929 | Frankenversand | München |
| 10893 | Königlich Essen | Brandenburg |
| 10891 | Lehmanns Marktstand | Frankfurt a.M. |
| 10878 | QUICK-Stop | Cunewalde |
| 10865 | QUICK-Stop | Cunewalde |
| 10862 | Lehmanns Marktstand | Frankfurt a.M. |
| 10859 | Frankenversand | München |
| 10853 | Blauer See Delikatessen | Mannheim |
| 10849 | Königlich Essen | Brandenburg |
| 10845 | QUICK-Stop | Cunewalde |
| 10835 | Alfred's Futterkiste | Berlin |
| 10833 | Ottilies Käseladen | Köln |
| 10825 | Drachenblut Delikatessen | Aachen |
| 10817 | Königlich Essen | Brandenburg |
| 10799 | Königlich Essen | Brandenburg |
| 10797 | Drachenblut Delikatessen | Aachen |
| 10791 | Frankenversand | München |
| 10788 | QUICK-Stop | Cunewalde |
| 10779 | Morgenstern Gesundkost | Leipzig |
| 10772 | Lehmanns Marktstand | Frankfurt a.M. |
| 10766 | Ottilies Käseladen | Köln |

| ID | Name | City |
|-------|-------------------------|----------------|
| 10765 | QUICK-Stop | Cunewalde |
| 10745 | QUICK-Stop | Cunewalde |
| 10721 | QUICK-Stop | Cunewalde |
| 10718 | Königlich Essen | Brandenburg |
| 10717 | Frankenversand | München |
| 10702 | Alfred's Futterkiste | Berlin |
| 10699 | Morgenstern Gesundkost | Leipzig |
| 10694 | QUICK-Stop | Cunewalde |
| 10692 | Alfred's Futterkiste | Berlin |
| 10691 | QUICK-Stop | Cunewalde |
| 10684 | Ottilies Käseladen | Köln |
| 10675 | Frankenversand | München |
| 10670 | Frankenversand | München |
| 10668 | Die Wandernde Kuh | Stuttgart |
| 10658 | QUICK-Stop | Cunewalde |
| 10653 | Frankenversand | München |
| 10651 | Die Wandernde Kuh | Stuttgart |
| 10643 | Alfreds Futterkiste | Berlin |
| 10640 | Die Wandernde Kuh | Stuttgart |
| 10632 | Die Wandernde Kuh | Stuttgart |
| 10630 | Königlich Essen | Brandenburg |
| 10623 | Frankenversand | München |
| 10614 | Blauer See Delikatessen | Mannheim |
| 10608 | Toms Spezialitäten | Münster |
| 10592 | Lehmanns Marktstand | Frankfurt a.M. |
| 10593 | Lehmanns Marktstand | Frankfurt a.M. |
| 10588 | QUICK-Stop | Cunewalde |
| 10582 | Blauer See Delikatessen | Mannheim |
| 10580 | Ottilies Käseladen | Köln |
| 10575 | Morgenstern Gesundkost | Leipzig |
| 10560 | Frankenversand | München |
| 10557 | Lehmanns Marktstand | Frankfurt a.M. |
| 10554 | Ottilies Käseladen | Köln |
| 10548 | Toms Spezialitäten | Münster |

| ID | Name | City |
|-------|--------------------------|----------------|
| 10549 | QUICK-Stop | Cunewalde |
| 10542 | Königlich Essen | Brandenburg |
| 10540 | QUICK-Stop | Cunewalde |
| 10536 | Lehmanns Marktstand | Frankfurt a.M. |
| 10534 | Lehmanns Marktstand | Frankfurt a.M. |
| 10527 | QUICK-Stop | Cunewalde |
| 10522 | Lehmanns Marktstand | Frankfurt a.M. |
| 10515 | QUICK-Stop | Cunewalde |
| 10513 | Die Wandernde Kuh | Stuttgart |
| 10508 | Ottilies Käseladen | Köln |
| 10509 | Blauer See Delikatessen | Mannheim |
| 10506 | Königlich Essen | Brandenburg |
| 10501 | Blauer See Delikatessen | Mannheim |
| 10497 | Lehmanns Marktstand | Frankfurt a.M. |
| 10488 | Frankenversand | München |
| 10468 | Königlich Essen | Brandenburg |
| 10456 | Königlich Essen | Brandenburg |
| 10457 | Königlich Essen | Brandenburg |
| 10451 | QUICK-Stop | Cunewalde |
| 10446 | Toms Spezialitäten | Münster |
| 10438 | Toms Spezialitäten | Münster |
| 10418 | QUICK-Stop | Cunewalde |
| 10407 | Ottilies Käseladen | Köln |
| 10396 | Frankenversand | München |
| 10391 | Drachenblut Delikatessen | Aachen |
| 10363 | Drachenblut Delikatessen | Aachen |
| 10361 | QUICK-Stop | Cunewalde |
| 10356 | Die Wandernde Kuh | Stuttgart |
| 10348 | Die Wandernde Kuh | Stuttgart |
| 10345 | QUICK-Stop | Cunewalde |
| 10342 | Frankenversand | München |
| 10343 | Lehmanns Marktstand | Frankfurt a.M. |
| 10337 | Frankenversand | München |
| 10325 | Königlich Essen | Brandenburg |

| ID | Name | City |
|-------|------------------------|----------------|
| 10323 | Königlich Essen | Brandenburg |
| 10312 | Die Wandernde Kuh | Stuttgart |
| 10313 | QUICK-Stop | Cunewalde |
| 10301 | Die Wandernde Kuh | Stuttgart |
| 10286 | QUICK-Stop | Cunewalde |
| 10284 | Lehmanns Marktstand | Frankfurt a.M. |
| 10285 | QUICK-Stop | Cunewalde |
| 10279 | Lehmanns Marktstand | Frankfurt a.M. |
| 10277 | Morgenstern Gesundkost | Leipzig |
| 10273 | QUICK-Stop | Cunewalde |
| 10267 | Frankenversand | München |
| 10260 | Ottilies Käseladen | Köln |
| 10249 | Toms Spezialitäten | Münster |

Count : 122

Countries Starting With Letter : I

10. Ireland

| | | |
|-------|------------------------------|------|
| 11063 | Hungry Owl All-Night Grocers | Cork |
| 10985 | Hungry Owl All-Night Grocers | Cork |
| 10912 | Hungry Owl All-Night Grocers | Cork |
| 10897 | Hungry Owl All-Night Grocers | Cork |
| 10736 | Hungry Owl All-Night Grocers | Cork |
| 10712 | Hungry Owl All-Night Grocers | Cork |
| 10701 | Hungry Owl All-Night Grocers | Cork |
| 10687 | Hungry Owl All-Night Grocers | Cork |
| 10661 | Hungry Owl All-Night Grocers | Cork |
| 10646 | Hungry Owl All-Night Grocers | Cork |
| 10567 | Hungry Owl All-Night Grocers | Cork |
| 10516 | Hungry Owl All-Night Grocers | Cork |
| 10503 | Hungry Owl All-Night Grocers | Cork |
| 10429 | Hungry Owl All-Night Grocers | Cork |
| 10380 | Hungry Owl All-Night Grocers | Cork |
| 10373 | Hungry Owl All-Night Grocers | Cork |
| 10335 | Hungry Owl All-Night Grocers | Cork |

| ID | Name | City |
|-------|------------------------------|------|
| 10309 | Hungry Owl All-Night Grocers | Cork |
| 10298 | Hungry Owl All-Night Grocers | Cork |

Count : 19

11. Italy

| | | |
|-------|------------------------------|---------------|
| 11062 | Reggiani Caseifici | Reggio Emilia |
| 11060 | Franchi S.p.A. | Torino |
| 11026 | Franchi S.p.A. | Torino |
| 11010 | Reggiani Caseifici | Reggio Emilia |
| 10950 | Magazzini Alimentari Riuniti | Bergamo |
| 10942 | Reggiani Caseifici | Reggio Emilia |
| 10939 | Magazzini Alimentari Riuniti | Bergamo |
| 10908 | Reggiani Caseifici | Reggio Emilia |
| 10818 | Magazzini Alimentari Riuniti | Bergamo |
| 10812 | Reggiani Caseifici | Reggio Emilia |
| 10807 | Franchi S.p.A. | Torino |
| 10784 | Magazzini Alimentari Riuniti | Bergamo |
| 10754 | Magazzini Alimentari Riuniti | Bergamo |
| 10753 | Franchi S.p.A. | Torino |
| 10727 | Reggiani Caseifici | Reggio Emilia |
| 10710 | Franchi S.p.A. | Torino |
| 10655 | Reggiani Caseifici | Reggio Emilia |
| 10635 | Magazzini Alimentari Riuniti | Bergamo |
| 10586 | Reggiani Caseifici | Reggio Emilia |
| 10562 | Reggiani Caseifici | Reggio Emilia |
| 10467 | Magazzini Alimentari Riuniti | Bergamo |
| 10443 | Reggiani Caseifici | Reggio Emilia |
| 10428 | Reggiani Caseifici | Reggio Emilia |
| 10422 | Franchi S.p.A. | Torino |
| 10404 | Magazzini Alimentari Riuniti | Bergamo |
| 10300 | Magazzini Alimentari Riuniti | Bergamo |
| 10288 | Reggiani Caseifici | Reggio Emilia |
| 10275 | Magazzini Alimentari Riuniti | Bergamo |

Count : 28

| ID | Name | City |
|---|------------------------------------|-------------------|
| <i>Countries Starting With Letter : M</i> | | |
| 12. Mexico | | |
| 11073 | Pericles Comidas clásicas | México D.F. |
| 11069 | Tortuga Restaurante | México D.F. |
| 10995 | Pericles Comidas clásicas | México D.F. |
| 10926 | Ana Trujillo Emparedados y helados | México D.F. |
| 10915 | Tortuga Restaurante | México D.F. |
| 10856 | Antonio Moreno Taquería | México D.F. |
| 10842 | Tortuga Restaurante | México D.F. |
| 10759 | Ana Trujillo Emparedados y helados | México D.F. |
| 10682 | Antonio Moreno Taquería | México D.F. |
| 10676 | Tortuga Restaurante | México D.F. |
| 10677 | Antonio Moreno Taquería | México D.F. |
| 10625 | Ana Trujillo Emparedados y helados | México D.F. |
| 10576 | Tortuga Restaurante | México D.F. |
| 10573 | Antonio Moreno Taquería | México D.F. |
| 10535 | Antonio Moreno Taquería | México D.F. |
| 10518 | Tortuga Restaurante | México D.F. |
| 10507 | Antonio Moreno Taquería | México D.F. |
| 10502 | Pericles Comidas clásicas | México D.F. |
| 10474 | Pericles Comidas clásicas | México D.F. |
| 10365 | Antonio Moreno Taquería | México D.F. |
| 10354 | Pericles Comidas clásicas | México D.F. |
| 10322 | Pericles Comidas clásicas | México D.F. |
| 10319 | Tortuga Restaurante | México D.F. |
| 10308 | Ana Trujillo Emparedados y helados | México D.F. |
| 10304 | Tortuga Restaurante | México D.F. |
| 10293 | Tortuga Restaurante | México D.F. |
| 10276 | Tortuga Restaurante | México D.F. |
| 10259 | Centro comercial Moctezuma | México D.F. |
| | | Count : 28 |

| ID | Name | City |
|----|------|------|
|----|------|------|

Countries Starting With Letter : N

13. Norway

| | | |
|-------|---------------|------------------|
| 11015 | Santé Gourmet | Stavern |
| 10909 | Santé Gourmet | Stavern |
| 10831 | Santé Gourmet | Stavern |
| 10639 | Santé Gourmet | Stavern |
| 10520 | Santé Gourmet | Stavern |
| 10387 | Santé Gourmet | Stavern |
| | | Count : 6 |

Countries Starting With Letter : P

14. Poland

| | | |
|-------|---------------|------------------|
| 11044 | Wolski Zajazd | Warszawa |
| 10998 | Wolski Zajazd | Warszawa |
| 10906 | Wolski Zajazd | Warszawa |
| 10870 | Wolski Zajazd | Warszawa |
| 10792 | Wolski Zajazd | Warszawa |
| 10611 | Wolski Zajazd | Warszawa |
| 10374 | Wolski Zajazd | Warszawa |
| | | Count : 7 |

15. Portugal

| | | |
|-------|--------------------------------|--------|
| 11007 | Princesa Isabel Vinhos | Lisboa |
| 10963 | Furia Bacalhau e Frutos do Mar | Lisboa |
| 10664 | Furia Bacalhau e Frutos do Mar | Lisboa |
| 10604 | Furia Bacalhau e Frutos do Mar | Lisboa |
| 10551 | Furia Bacalhau e Frutos do Mar | Lisboa |
| 10491 | Furia Bacalhau e Frutos do Mar | Lisboa |
| 10477 | Princesa Isabel Vinhos | Lisboa |
| 10464 | Furia Bacalhau e Frutos do Mar | Lisboa |
| 10433 | Princesa Isabel Vinhos | Lisboa |
| 10397 | Princesa Isabel Vinhos | Lisboa |
| 10352 | Furia Bacalhau e Frutos do Mar | Lisboa |
| 10336 | Princesa Isabel Vinhos | Lisboa |
| 10328 | Furia Bacalhau e Frutos do Mar | Lisboa |

| ID | Name | City |
|----|------|------|
|----|------|------|

Count : 13

Countries Starting With Letter : S

16. Spain

| | | |
|-------|---------------------------|-----------|
| 11037 | Godos Cocina Típica | Sevilla |
| 11013 | Romero y tomillo | Madrid |
| 11009 | Godos Cocina Típica | Sevilla |
| 10970 | Bólido Comidas preparadas | Madrid |
| 10948 | Godos Cocina Típica | Sevilla |
| 10928 | Galería del gastronómo | Barcelona |
| 10917 | Romero y tomillo | Madrid |
| 10911 | Godos Cocina Típica | Sevilla |
| 10888 | Godos Cocina Típica | Sevilla |
| 10887 | Galería del gastronómo | Barcelona |
| 10874 | Godos Cocina Típica | Sevilla |
| 10872 | Godos Cocina Típica | Sevilla |
| 10801 | Bólido Comidas preparadas | Madrid |
| 10629 | Godos Cocina Típica | Sevilla |
| 10568 | Galería del gastronómo | Barcelona |
| 10550 | Godos Cocina Típica | Sevilla |
| 10426 | Galería del gastronómo | Barcelona |
| 10366 | Galería del gastronómo | Barcelona |
| 10326 | Bólido Comidas preparadas | Madrid |
| 10306 | Romero y tomillo | Madrid |
| 10303 | Godos Cocina Típica | Sevilla |
| 10282 | Romero y tomillo | Madrid |
| 10281 | Romero y tomillo | Madrid |

Count : 23

17. Sweden

| | | |
|-------|----------------|--------|
| 11050 | Folk och fä HB | Bräcke |
| 11001 | Folk och fä HB | Bräcke |
| 10993 | Folk och fä HB | Bräcke |
| 10980 | Folk och fä HB | Bräcke |
| 10977 | Folk och fä HB | Bräcke |

| ID | Name | City |
|-------|--------------------|--------|
| 10955 | Folk och fä HB | Bräcke |
| 10924 | Berglunds snabbköp | Luleå |
| 10902 | Folk och fä HB | Bräcke |
| 10880 | Folk och fä HB | Bräcke |
| 10875 | Berglunds snabbköp | Luleå |
| 10866 | Berglunds snabbköp | Luleå |
| 10857 | Berglunds snabbköp | Luleå |
| 10837 | Berglunds snabbköp | Luleå |
| 10824 | Folk och fä HB | Bräcke |
| 10778 | Berglunds snabbköp | Luleå |
| 10774 | Folk och fä HB | Bräcke |
| 10762 | Folk och fä HB | Bräcke |
| 10733 | Berglunds snabbköp | Luleå |
| 10703 | Folk och fä HB | Bräcke |
| 10689 | Berglunds snabbköp | Luleå |
| 10672 | Berglunds snabbköp | Luleå |
| 10654 | Berglunds snabbköp | Luleå |
| 10626 | Berglunds snabbköp | Luleå |
| 10572 | Berglunds snabbköp | Luleå |
| 10561 | Folk och fä HB | Bräcke |
| 10533 | Folk och fä HB | Bräcke |
| 10524 | Berglunds snabbköp | Luleå |
| 10460 | Folk och fä HB | Bräcke |
| 10444 | Berglunds snabbköp | Luleå |
| 10445 | Berglunds snabbköp | Luleå |
| 10434 | Folk och fä HB | Bräcke |
| 10384 | Berglunds snabbköp | Luleå |
| 10378 | Folk och fä HB | Bräcke |
| 10327 | Folk och fä HB | Bräcke |
| 10280 | Berglunds snabbköp | Luleå |
| 10278 | Berglunds snabbköp | Luleå |
| 10264 | Folk och fä HB | Bräcke |

Count : 37

| ID | Name | City |
|------------------------|--------------------|-------------------|
| 18. Switzerland | | |
| 11075 | Richter Supermarkt | Genève |
| 11041 | Chop-suey Chinese | Bern |
| 11033 | Richter Supermarkt | Genève |
| 11029 | Chop-suey Chinese | Bern |
| 10966 | Chop-suey Chinese | Bern |
| 10951 | Richter Supermarkt | Genève |
| 10931 | Richter Supermarkt | Genève |
| 10758 | Richter Supermarkt | Genève |
| 10751 | Richter Supermarkt | Genève |
| 10746 | Chop-suey Chinese | Bern |
| 10731 | Chop-suey Chinese | Bern |
| 10666 | Richter Supermarkt | Genève |
| 10537 | Richter Supermarkt | Genève |
| 10519 | Chop-suey Chinese | Bern |
| 10419 | Richter Supermarkt | Genève |
| 10370 | Chop-suey Chinese | Bern |
| 10254 | Chop-suey Chinese | Bern |
| 10255 | Richter Supermarkt | Genève |
| | | Count : 18 |

Countries Starting With Letter : U

| | | |
|---------------|--------------------|------------|
| 19. UK | | |
| 11056 | Eastern Connection | London |
| 11057 | North/South | London |
| 11047 | Eastern Connection | London |
| 11024 | Eastern Connection | London |
| 11023 | B's Beverages | London |
| 11016 | Around the Horn | Colchester |
| 10987 | Eastern Connection | London |
| 10953 | Around the Horn | Colchester |
| 10947 | B's Beverages | London |
| 10943 | B's Beverages | London |
| 10933 | Island Trading | Cowes |

| ID | Name | City |
|-------|-----------------------|------------|
| 10920 | Around the Horn | Colchester |
| 10869 | Seven Seas Imports | London |
| 10864 | Around the Horn | Colchester |
| 10848 | Consolidated Holdings | London |
| 10829 | Island Trading | Cowes |
| 10804 | Seven Seas Imports | London |
| 10800 | Seven Seas Imports | London |
| 10798 | Island Trading | Cowes |
| 10793 | Around the Horn | Colchester |
| 10768 | Around the Horn | Colchester |
| 10752 | North/South | London |
| 10749 | Island Trading | Cowes |
| 10743 | Around the Horn | Colchester |
| 10741 | Around the Horn | Colchester |
| 10726 | Eastern Connection | London |
| 10707 | Around the Horn | Colchester |
| 10674 | Island Trading | Cowes |
| 10621 | Island Trading | Cowes |
| 10599 | B's Beverages | London |
| 10578 | B's Beverages | London |
| 10558 | Around the Horn | Colchester |
| 10547 | Seven Seas Imports | London |
| 10538 | B's Beverages | London |
| 10539 | B's Beverages | London |
| 10532 | Eastern Connection | London |
| 10523 | Seven Seas Imports | London |
| 10517 | North/South | London |
| 10484 | B's Beverages | London |
| 10472 | Seven Seas Imports | London |
| 10473 | Island Trading | Cowes |
| 10471 | B's Beverages | London |
| 10462 | Consolidated Holdings | London |
| 10453 | Around the Horn | Colchester |
| 10435 | Consolidated Holdings | London |

| ID | Name | City |
|-------|--------------------|------------|
| 10400 | Eastern Connection | London |
| 10388 | Seven Seas Imports | London |
| 10383 | Around the Horn | Colchester |
| 10377 | Seven Seas Imports | London |
| 10364 | Eastern Connection | London |
| 10359 | Seven Seas Imports | London |
| 10355 | Around the Horn | Colchester |
| 10321 | Island Trading | Cowes |
| 10318 | Island Trading | Cowes |
| 10315 | Island Trading | Cowes |
| 10289 | B's Beverages | London |

Count : 56

20. USA

| | | |
|-------|----------------------------|-------------|
| 11077 | Rattlesnake Canyon Grocery | Albuquerque |
| 11066 | White Clover Markets | Seattle |
| 11064 | Save-a-lot Markets | Boise |
| 11061 | Great Lakes Food Market | Eugene |
| 11040 | Great Lakes Food Market | Eugene |
| 11034 | Old World Delicatessen | Anchorage |
| 11032 | White Clover Markets | Seattle |
| 11030 | Save-a-lot Markets | Boise |
| 11031 | Save-a-lot Markets | Boise |
| 11018 | Lonesome Pine Restaurant | Portland |
| 11006 | Great Lakes Food Market | Eugene |
| 11002 | Save-a-lot Markets | Boise |
| 11003 | The Cracker Box | Butte |
| 11000 | Rattlesnake Canyon Grocery | Albuquerque |
| 10992 | The Big Cheese | Portland |
| 10988 | Rattlesnake Canyon Grocery | Albuquerque |
| 10984 | Save-a-lot Markets | Boise |
| 10983 | Save-a-lot Markets | Boise |
| 10974 | Split Rail Beer & Ale | Lander |
| 10965 | Old World Delicatessen | Anchorage |
| 10941 | Save-a-lot Markets | Boise |

| ID | Name | City |
|-------|-----------------------------------|---------------|
| 10936 | Great Lakes Food Market | Eugene |
| 10904 | White Clover Markets | Seattle |
| 10894 | Save-a-lot Markets | Boise |
| 10889 | Rattlesnake Canyon Grocery | Albuquerque |
| 10884 | Let's Stop N Shop | San Francisco |
| 10882 | Save-a-lot Markets | Boise |
| 10883 | Lonesome Pine Restaurant | Portland |
| 10867 | Lonesome Pine Restaurant | Portland |
| 10861 | White Clover Markets | Seattle |
| 10855 | Old World Delicatessen | Anchorage |
| 10852 | Rattlesnake Canyon Grocery | Albuquerque |
| 10847 | Save-a-lot Markets | Boise |
| 10822 | Trail's Head Gourmet Provisioners | Kirkland |
| 10820 | Rattlesnake Canyon Grocery | Albuquerque |
| 10821 | Split Rail Beer & Ale | Lander |
| 10816 | Great Lakes Food Market | Eugene |
| 10815 | Save-a-lot Markets | Boise |
| 10808 | Old World Delicatessen | Anchorage |
| 10805 | The Big Cheese | Portland |
| 10775 | The Cracker Box | Butte |
| 10761 | Rattlesnake Canyon Grocery | Albuquerque |
| 10756 | Split Rail Beer & Ale | Lander |
| 10757 | Save-a-lot Markets | Boise |
| 10748 | Save-a-lot Markets | Boise |
| 10740 | White Clover Markets | Seattle |
| 10735 | Let's Stop N Shop | San Francisco |
| 10722 | Save-a-lot Markets | Boise |
| 10723 | White Clover Markets | Seattle |
| 10719 | Let's Stop N Shop | San Francisco |
| 10714 | Save-a-lot Markets | Boise |
| 10713 | Save-a-lot Markets | Boise |
| 10711 | Save-a-lot Markets | Boise |
| 10708 | The Big Cheese | Portland |
| 10706 | Old World Delicatessen | Anchorage |

| ID | Name | City |
|-------|-----------------------------------|---------------|
| 10700 | Save-a-lot Markets | Boise |
| 10696 | White Clover Markets | Seattle |
| 10693 | White Clover Markets | Seattle |
| 10680 | Old World Delicatessen | Anchorage |
| 10681 | Great Lakes Food Market | Eugene |
| 10678 | Save-a-lot Markets | Boise |
| 10665 | Lonesome Pine Restaurant | Portland |
| 10662 | Lonesome Pine Restaurant | Portland |
| 10660 | Hungry Coyote Import Store | Elgin |
| 10656 | Great Lakes Food Market | Eugene |
| 10657 | Save-a-lot Markets | Boise |
| 10627 | Save-a-lot Markets | Boise |
| 10624 | The Cracker Box | Butte |
| 10616 | Great Lakes Food Market | Eugene |
| 10617 | Great Lakes Food Market | Eugene |
| 10612 | Save-a-lot Markets | Boise |
| 10607 | Save-a-lot Markets | Boise |
| 10603 | Save-a-lot Markets | Boise |
| 10600 | Hungry Coyote Import Store | Elgin |
| 10598 | Rattlesnake Canyon Grocery | Albuquerque |
| 10596 | White Clover Markets | Seattle |
| 10594 | Old World Delicatessen | Anchorage |
| 10589 | Great Lakes Food Market | Eugene |
| 10579 | Let's Stop N Shop | San Francisco |
| 10577 | Trail's Head Gourmet Provisioners | Kirkland |
| 10574 | Trail's Head Gourmet Provisioners | Kirkland |
| 10569 | Rattlesnake Canyon Grocery | Albuquerque |
| 10564 | Rattlesnake Canyon Grocery | Albuquerque |
| 10555 | Save-a-lot Markets | Boise |
| 10544 | Lonesome Pine Restaurant | Portland |
| 10545 | Lazy K Kountry Store | Walla Walla |
| 10528 | Great Lakes Food Market | Eugene |
| 10510 | Save-a-lot Markets | Boise |
| 10504 | White Clover Markets | Seattle |

| ID | Name | City |
|-------|----------------------------|-------------|
| 10482 | Lazy K Kountry Store | Walla Walla |
| 10483 | White Clover Markets | Seattle |
| 10479 | Rattlesnake Canyon Grocery | Albuquerque |
| 10469 | White Clover Markets | Seattle |
| 10452 | Save-a-lot Markets | Boise |
| 10440 | Save-a-lot Markets | Boise |
| 10441 | Old World Delicatessen | Anchorage |
| 10432 | Split Rail Beer & Ale | Lander |
| 10415 | Hungry Coyote Import Store | Elgin |
| 10401 | Rattlesnake Canyon Grocery | Albuquerque |
| 10398 | Save-a-lot Markets | Boise |
| 10394 | Hungry Coyote Import Store | Elgin |
| 10393 | Save-a-lot Markets | Boise |
| 10385 | Split Rail Beer & Ale | Lander |
| 10375 | Hungry Coyote Import Store | Elgin |
| 10369 | Split Rail Beer & Ale | Lander |
| 10349 | Split Rail Beer & Ale | Lander |
| 10346 | Rattlesnake Canyon Grocery | Albuquerque |
| 10344 | White Clover Markets | Seattle |
| 10338 | Old World Delicatessen | Anchorage |
| 10329 | Split Rail Beer & Ale | Lander |
| 10324 | Save-a-lot Markets | Boise |
| 10316 | Rattlesnake Canyon Grocery | Albuquerque |
| 10317 | Lonesome Pine Restaurant | Portland |
| 10314 | Rattlesnake Canyon Grocery | Albuquerque |
| 10310 | The Big Cheese | Portland |
| 10307 | Lonesome Pine Restaurant | Portland |
| 10305 | Old World Delicatessen | Anchorage |
| 10294 | Rattlesnake Canyon Grocery | Albuquerque |
| 10272 | Rattlesnake Canyon Grocery | Albuquerque |
| 10271 | Split Rail Beer & Ale | Lander |
| 10269 | White Clover Markets | Seattle |
| 10262 | Rattlesnake Canyon Grocery | Albuquerque |

Count : 122

| ID | Name | City |
|---|----------------------|-----------------|
| <i>Countries Starting With Letter : V</i> | | |
| 21. Venezuela | | |
| 11071 | LILA-Supermercado | Barquisimeto |
| 11065 | LILA-Supermercado | Barquisimeto |
| 11055 | HILARION-Abastos | San Cristóbal |
| 11039 | LINO-Delicateses | I. de Margarita |
| 11014 | LINO-Delicateses | I. de Margarita |
| 10997 | LILA-Supermercado | Barquisimeto |
| 10976 | HILARION-Abastos | San Cristóbal |
| 10960 | HILARION-Abastos | San Cristóbal |
| 10957 | HILARION-Abastos | San Cristóbal |
| 10954 | LINO-Delicateses | I. de Margarita |
| 10919 | LINO-Delicateses | I. de Margarita |
| 10901 | HILARION-Abastos | San Cristóbal |
| 10899 | LILA-Supermercado | Barquisimeto |
| 10863 | HILARION-Abastos | San Cristóbal |
| 10840 | LINO-Delicateses | I. de Margarita |
| 10838 | LINO-Delicateses | I. de Margarita |
| 10823 | LILA-Supermercado | Barquisimeto |
| 10811 | LINO-Delicateses | I. de Margarita |
| 10796 | HILARION-Abastos | San Cristóbal |
| 10785 | GROSELLA-Restaurante | Caracas |
| 10780 | LILA-Supermercado | Barquisimeto |
| 10729 | LINO-Delicateses | I. de Margarita |
| 10705 | HILARION-Abastos | San Cristóbal |
| 10697 | LINO-Delicateses | I. de Margarita |
| 10641 | HILARION-Abastos | San Cristóbal |
| 10638 | LINO-Delicateses | I. de Margarita |
| 10613 | HILARION-Abastos | San Cristóbal |
| 10601 | HILARION-Abastos | San Cristóbal |
| 10552 | HILARION-Abastos | San Cristóbal |
| 10543 | LILA-Supermercado | Barquisimeto |
| 10498 | HILARION-Abastos | San Cristóbal |

| ID | Name | City |
|-------|----------------------|-----------------|
| 10499 | LILA-Supermercado | Barquisimeto |
| 10490 | HILARION-Abastos | San Cristóbal |
| 10486 | HILARION-Abastos | San Cristóbal |
| 10485 | LINO-Delicateses | I. de Margarita |
| 10476 | HILARION-Abastos | San Cristóbal |
| 10461 | LILA-Supermercado | Barquisimeto |
| 10405 | LINO-Delicateses | I. de Margarita |
| 10395 | HILARION-Abastos | San Cristóbal |
| 10381 | LILA-Supermercado | Barquisimeto |
| 10357 | LILA-Supermercado | Barquisimeto |
| 10330 | LILA-Supermercado | Barquisimeto |
| 10296 | LILA-Supermercado | Barquisimeto |
| 10283 | LILA-Supermercado | Barquisimeto |
| 10268 | GROSELLA-Restaurante | Caracas |
| 10257 | HILARION-Abastos | San Cristóbal |

Count : 46

There were 830 orders on this report.