

There are 253 orders on  
this report, with a total  
freight of 16410.71

## Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **A**

**Argentina**

**Mon, Feb 17, 1997**

<b>Order</b>	10448	<b>Order</b>	10409
<b>Name</b>	Rancho grande	<b>Name</b>	Océano Atlántico Ltda.
<b>City</b>	Buenos Aires	<b>City</b>	Buenos Aires
<b>Region</b>		<b>Region</b>	
<b>Date</b>	17/02/1997	<b>Date</b>	09/01/1997
<b>Freight</b>	38.82	<b>Freight</b>	29.83

**Count : 2 Total : 68.65**

**Austria**

**Fri, Mar 28, 1997**

<b>Order</b>	10489	<b>Order</b>	10442	<b>Order</b>	<b>10430</b>
<b>Name</b>	Piccolo und mehr	<b>Name</b>	Ernst Handel	<b>Name</b>	Ernst Handel
<b>City</b>	Salzburg	<b>City</b>	Graz	<b>City</b>	Graz
<b>Region</b>		<b>Region</b>		<b>Region</b>	
<b>Date</b>	28/03/1997	<b>Date</b>	11/02/1997	<b>Date</b>	30/01/1997
<b>Freight</b>	5.29	<b>Freight</b>	47.94	<b>Freight</b>	458.78
<b>Order</b>	10427	<b>Order</b>	10403	<b>Order</b>	10402
<b>Name</b>	Piccolo und mehr	<b>Name</b>	Ernst Handel	<b>Name</b>	Ernst Handel
<b>City</b>	Salzburg	<b>City</b>	Graz	<b>City</b>	Graz
<b>Region</b>		<b>Region</b>		<b>Region</b>	
<b>Date</b>	27/01/1997	<b>Date</b>	03/01/1997	<b>Date</b>	02/01/1997
<b>Freight</b>	31.29	<b>Freight</b>	73.79	<b>Freight</b>	67.88
<b>Order</b>	10392	<b>Order</b>	<b>10390</b>	<b>Order</b>	10382
<b>Name</b>	Piccolo und mehr	<b>Name</b>	Ernst Handel	<b>Name</b>	Ernst Handel
<b>City</b>	Salzburg	<b>City</b>	Graz	<b>City</b>	Graz
<b>Region</b>		<b>Region</b>		<b>Region</b>	
<b>Date</b>	24/12/1996	<b>Date</b>	23/12/1996	<b>Date</b>	13/12/1996
<b>Freight</b>	122.46	<b>Freight</b>	126.38	<b>Freight</b>	94.77
<b>Order</b>	10368	<b>Order</b>	10353	<b>Order</b>	10351
<b>Name</b>	Ernst Handel	<b>Name</b>	Piccolo und mehr	<b>Name</b>	Ernst Handel
<b>City</b>	Graz	<b>City</b>	Salzburg	<b>City</b>	Graz
<b>Region</b>		<b>Region</b>		<b>Region</b>	
<b>Date</b>	29/11/1996	<b>Date</b>	13/11/1996	<b>Date</b>	11/11/1996
<b>Freight</b>	101.95	<b>Freight</b>	360.63	<b>Freight</b>	162.33
<b>Order</b>	10263	<b>Order</b>	10258		
<b>Name</b>	Ernst Handel	<b>Name</b>	Ernst Handel		
<b>City</b>	Graz	<b>City</b>	Graz		
<b>Region</b>		<b>Region</b>			
<b>Date</b>	23/07/1996	<b>Date</b>	17/07/1996		
<b>Freight</b>	146.06	<b>Freight</b>	140.51		

**Count : 16 Total : 2008.71**

**Count : 16 Total : 2008.71**

**Count : 16 Total : 2008.71**

**Count : 16**

**1/20**

**Total : 2008.71**

Northwind Order List

Max order ID is : 10500

Orders

Orders

Orders

Countries Starting With Letter : A

Austria		Fri, Mar 28, 1997	
Count :	14	Total :	1940.06
Count :	16	Total :	2008.71

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **B**

**Belgium**

**Fri, Mar 14, 1997**

<b>Order</b> 10475 <b>Name</b> Suprêmes délices <b>City</b> Charleroi <b>Region</b> <b>Date</b> 14/03/1997 <b>Freight</b> 68.52	<b>Order</b> 10463 <b>Name</b> Suprêmes délices <b>City</b> Charleroi <b>Region</b> <b>Date</b> 04/03/1997 <b>Freight</b> 14.78	<b>Order</b> 10458 <b>Name</b> Suprêmes délices <b>City</b> Charleroi <b>Region</b> <b>Date</b> 26/02/1997 <b>Freight</b> 147.06
<b>Order</b> 10302 <b>Name</b> Suprêmes délices <b>City</b> Charleroi <b>Region</b> <b>Date</b> 10/09/1996 <b>Freight</b> 6.27	<b>Order</b> 10252 <b>Name</b> Suprêmes délices <b>City</b> Charleroi <b>Region</b> <b>Date</b> 09/07/1996 <b>Freight</b> 51.30	
<b>Count :</b> 5 <b>Total :</b> 287.93		

**Brazil**

**Fri, Apr 4, 1997**

<b>Order</b> 10496 <b>Name</b> Tradição Hipermercados <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 04/04/1997 <b>Freight</b> 46.77	<b>Order</b> 10494 <b>Name</b> Comércio Mineiro <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 02/04/1997 <b>Freight</b> 65.99	<b>Order</b> 10487 <b>Name</b> Queen Cozinha <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 26/03/1997 <b>Freight</b> 71.07
<b>Order</b> 10481 <b>Name</b> Ricardo Adocicados <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 20/03/1997 <b>Freight</b> 64.33	<b>Order</b> 10466 <b>Name</b> Comércio Mineiro <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 06/03/1997 <b>Freight</b> 11.93	<b>Order</b> 10447 <b>Name</b> Ricardo Adocicados <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 14/02/1997 <b>Freight</b> 68.66
<b>Order</b> 10423 <b>Name</b> Gourmet Lanchonetes <b>City</b> Campinas <b>Region</b> SP <b>Date</b> 23/01/1997 <b>Freight</b> 24.50	<b>Order</b> 10421 <b>Name</b> Que Delícia <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 21/01/1997 <b>Freight</b> 99.23	<b>Order</b> 10420 <b>Name</b> Wellington Importadora <b>City</b> Resende <b>Region</b> SP <b>Date</b> 21/01/1997 <b>Freight</b> 44.12
<b>Order</b> 10414 <b>Name</b> Família Arquibaldo <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 14/01/1997 <b>Freight</b> 21.48	<b>Order</b> 10406 <b>Name</b> Queen Cozinha <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 07/01/1997 <b>Freight</b> 108.04	<b>Order</b> 10386 <b>Name</b> Família Arquibaldo <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 18/12/1996 <b>Freight</b> 13.99
<b>Order</b> 10379 <b>Name</b> Que Delícia <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 11/12/1996 <b>Freight</b> 45.03	<b>Order</b> 10372 <b>Name</b> Queen Cozinha <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 04/12/1996 <b>Freight</b> 890.78	<b>Order</b> 10347 <b>Name</b> Família Arquibaldo <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 06/11/1996 <b>Freight</b> 3.10

<b>Count :</b> 20 <b>Total :</b> 1866.95	<b>Count :</b> 20 <b>Total :</b> 1866.95	<b>Count :</b> 20 <b>Total :</b> 1866.95
--	--	--

**Count :** 20

**3/20**

**Total :** 1866.95

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **B**

**Brazil**

**Fri, Apr 4, 1997**

<b>Order</b> 10299 <b>Name</b> Ricardo Adocicados <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 06/09/1996 <b>Freight</b> 29.76	<b>Order</b> 10292 <b>Name</b> Tradiçao Hipermercados <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 28/08/1996 <b>Freight</b> 1.35	<b>Order</b> 10291 <b>Name</b> Que Delícia <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 27/08/1996 <b>Freight</b> 6.40
<b>Order</b> <b>10290</b> <b>Name</b> Comércio Mineiro <b>City</b> Sao Paulo <b>Region</b> SP <b>Date</b> 27/08/1996 <b>Freight</b> 79.70	<b>Order</b> 10287 <b>Name</b> Ricardo Adocicados <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 22/08/1996 <b>Freight</b> 12.76	<b>Order</b> 10261 <b>Name</b> Que Delícia <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 19/07/1996 <b>Freight</b> 3.05
<b>Order</b> 10256 <b>Name</b> Wellington Importadora <b>City</b> Resende <b>Region</b> SP <b>Date</b> 15/07/1996 <b>Freight</b> 13.97	<b>Order</b> 10253 <b>Name</b> Hanari Carnes <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 10/07/1996 <b>Freight</b> 58.17	<b>Order</b> <b>10250</b> <b>Name</b> Hanari Carnes <b>City</b> Rio de Janeiro <b>Region</b> RJ <b>Date</b> 08/07/1996 <b>Freight</b> 65.83

**Count : 24      Total : 1850.01**

**Count : 29      Total : 2137.94**

**Count : 9      Total : 270.99**

**Count : 9      Total : 270.99**

**Count : 9      Total : 270.99**

**Count : 9      4/20      Total : 270.99**

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **C**

Canada

Thu, Apr 3, 1997

<b>Order</b> 10495 <b>Name</b> Laughing Bacchus Wine Cellars <b>City</b> Vancouver <b>Region</b> BC <b>Date</b> 03/04/1997 <b>Freight</b> 4.65	<b>Order</b> 10492 <b>Name</b> Bottom-Dollar Markets <b>City</b> Tsawassen <b>Region</b> BC <b>Date</b> 01/04/1997 <b>Freight</b> 62.89	<b>Order</b> 10439 <b>Name</b> Mère Paillarde <b>City</b> Montréal <b>Region</b> Québec <b>Date</b> 07/02/1997 <b>Freight</b> 4.07
<b>Order</b> 10431 <b>Name</b> Bottom-Dollar Markets <b>City</b> Tsawassen <b>Region</b> BC <b>Date</b> 30/01/1997 <b>Freight</b> 44.17	<b>Order</b> 10424 <b>Name</b> Mère Paillarde <b>City</b> Montréal <b>Region</b> Québec <b>Date</b> 23/01/1997 <b>Freight</b> 370.61	<b>Order</b> 10411 <b>Name</b> Bottom-Dollar Markets <b>City</b> Tsawassen <b>Region</b> BC <b>Date</b> 10/01/1997 <b>Freight</b> 23.65
<b>Order</b> <b>10410</b> <b>Name</b> Bottom-Dollar Markets <b>City</b> Tsawassen <b>Region</b> BC <b>Date</b> 10/01/1997 <b>Freight</b> 2.40	<b>Order</b> 10389 <b>Name</b> Bottom-Dollar Markets <b>City</b> Tsawassen <b>Region</b> BC <b>Date</b> 20/12/1996 <b>Freight</b> 47.42	<b>Order</b> 10376 <b>Name</b> Mère Paillarde <b>City</b> Montréal <b>Region</b> Québec <b>Date</b> 09/12/1996 <b>Freight</b> 20.39
<b>Order</b> 10339 <b>Name</b> Mère Paillarde <b>City</b> Montréal <b>Region</b> Québec <b>Date</b> 28/10/1996 <b>Freight</b> 15.66	<b>Order</b> 10332 <b>Name</b> Mère Paillarde <b>City</b> Montréal <b>Region</b> Québec <b>Date</b> 17/10/1996 <b>Freight</b> 52.84	

Count : 11 Total : 648.75

Count : 11 Total : 648.75

Count : 11 Total : 648.75

Count : 11 Total : 648.75

Count : 11 Total : 648.75

Count : 11

5/20

Total : 648.75

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **D**

**Denmark**

**Wed, Mar 5, 1997**

<b>Order</b> 10465 <b>Name</b> Vaffeljernet <b>City</b> Århus <b>Region</b> <b>Date</b> 05/03/1997 <b>Freight</b> 145.04	<b>Order</b> 10417 <b>Name</b> Simons bistro <b>City</b> Kobenhavn <b>Region</b> <b>Date</b> 16/01/1997 <b>Freight</b> 70.29	<b>Order</b> 10399 <b>Name</b> Vaffeljernet <b>City</b> Århus <b>Region</b> <b>Date</b> 31/12/1996 <b>Freight</b> 27.36
<b>Order</b> 10367 <b>Name</b> Vaffeljernet <b>City</b> Århus <b>Region</b> <b>Date</b> 28/11/1996 <b>Freight</b> 13.55	<b>Order</b> 10341 <b>Name</b> Simons bistro <b>City</b> Kobenhavn <b>Region</b> <b>Date</b> 29/10/1996 <b>Freight</b> 26.78	

**Count :** 5 **Total :** 283.02

**Count :** 5 **Total :** 283.02

**Count :** 5 **Total :** 283.02

**Count :** 5 **Total :** 283.02

**Count :** 5 **Total :** 283.02

**Count :** 5 **6/20** **Total :** 283.02

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **F**

**Finland**

**Mon, Feb 24, 1997**

<b>Order</b> 10455 <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 24/02/1997 <b>Freight</b> 180.45	<b>Order</b> 10437 <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 05/02/1997 <b>Freight</b> 19.97	<b>Order</b> 10416 <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 16/01/1997 <b>Freight</b> 22.72
<b>Order</b> 10412 <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 13/01/1997 <b>Freight</b> 3.77	<b>Order</b> 10333 <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 18/10/1996 <b>Freight</b> 0.59	<b>Order</b> <b>10320</b> <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 03/10/1996 <b>Freight</b> 34.57
<b>Order</b> <b>10270</b> <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 01/08/1996 <b>Freight</b> 136.54	<b>Order</b> 10266 <b>Name</b> Wartian Herkku <b>City</b> Oulu <b>Region</b> <b>Date</b> 26/07/1996 <b>Freight</b> 25.73	

**Count :** 8 **Total :** 424.34

**France**

**Wed, Apr 9, 1997**

<b>Order</b> <b>10500</b> <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 09/04/1997 <b>Freight</b> 42.68	<b>Order</b> 10493 <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 02/04/1997 <b>Freight</b> 10.64	<b>Order</b> <b>10480</b> <b>Name</b> Folies gourmandes <b>City</b> Lille <b>Region</b> <b>Date</b> 20/03/1997 <b>Freight</b> 1.35
<b>Order</b> 10478 <b>Name</b> Victuailles en stock <b>City</b> Lyon <b>Region</b> <b>Date</b> 18/03/1997 <b>Freight</b> 4.81	<b>Order</b> <b>10470</b> <b>Name</b> Bon app' <b>City</b> Marseille <b>Region</b> <b>Date</b> 11/03/1997 <b>Freight</b> 64.56	<b>Order</b> 10459 <b>Name</b> Victuailles en stock <b>City</b> Lyon <b>Region</b> <b>Date</b> 27/02/1997 <b>Freight</b> 25.09
<b>Order</b> 10454 <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 21/02/1997 <b>Freight</b> 2.74	<b>Order</b> <b>10450</b> <b>Name</b> Victuailles en stock <b>City</b> Lyon <b>Region</b> <b>Date</b> 19/02/1997 <b>Freight</b> 7.23	<b>Order</b> 10449 <b>Name</b> Blondel père et fils <b>City</b> Strasbourg <b>Region</b> <b>Date</b> 18/02/1997 <b>Freight</b> 53.30
<b>Order</b> 10436 <b>Name</b> Blondel père et fils <b>City</b> Strasbourg <b>Region</b> <b>Date</b> 05/02/1997 <b>Freight</b> 156.66	<b>Order</b> 10425 <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 24/01/1997 <b>Freight</b> 7.93	<b>Order</b> 10413 <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 14/01/1997 <b>Freight</b> 95.66

**Count :** 20 **Total :** 896.99

**Count :** 20 **Total :** 896.99

**Count :** 20 **Total :** 896.99

**Count :** 20

**7/20**

**Total :** 896.99

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **F**

France

Wed, Apr 9, 1997

<b>Order</b> 10408 <b>Name</b> Folies gourmandes <b>City</b> Lille <b>Region</b> <b>Date</b> 08/01/1997 <b>Freight</b> 11.26	<b>Order</b> 10371 <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 03/12/1996 <b>Freight</b> 0.45	<b>Order</b> 10362 <b>Name</b> Bon app' <b>City</b> Marseille <b>Region</b> <b>Date</b> 25/11/1996 <b>Freight</b> 96.04
<b>Order</b> <b>10360</b> <b>Name</b> Blondel père et fils <b>City</b> Strasbourg <b>Region</b> <b>Date</b> 22/11/1996 <b>Freight</b> 131.70	<b>Order</b> 10358 <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 20/11/1996 <b>Freight</b> 19.64	<b>Order</b> <b>10350</b> <b>Name</b> La maison d'Asie <b>City</b> Toulouse <b>Region</b> <b>Date</b> 11/11/1996 <b>Freight</b> 64.19
<b>Order</b> <b>10340</b> <b>Name</b> Bon app' <b>City</b> Marseille <b>Region</b> <b>Date</b> 29/10/1996 <b>Freight</b> 166.31	<b>Order</b> 10334 <b>Name</b> Victuailles en stock <b>City</b> Lyon <b>Region</b> <b>Date</b> 21/10/1996 <b>Freight</b> 8.56	<b>Order</b> 10331 <b>Name</b> Bon app' <b>City</b> Marseille <b>Region</b> <b>Date</b> 16/10/1996 <b>Freight</b> 10.19
<b>Order</b> 10311 <b>Name</b> Du monde entier <b>City</b> Nantes <b>Region</b> <b>Date</b> 20/09/1996 <b>Freight</b> 24.69	<b>Order</b> 10297 <b>Name</b> Blondel père et fils <b>City</b> Strasbourg <b>Region</b> <b>Date</b> 04/09/1996 <b>Freight</b> 5.74	<b>Order</b> 10295 <b>Name</b> Vins et alcools Chevalier <b>City</b> Reims <b>Region</b> <b>Date</b> 02/09/1996 <b>Freight</b> 1.15
<b>Order</b> 10274 <b>Name</b> Vins et alcools Chevalier <b>City</b> Reims <b>Region</b> <b>Date</b> 06/08/1996 <b>Freight</b> 6.01	<b>Order</b> 10265 <b>Name</b> Blondel père et fils <b>City</b> Strasbourg <b>Region</b> <b>Date</b> 25/07/1996 <b>Freight</b> 55.28	<b>Order</b> 10251 <b>Name</b> Victuailles en stock <b>City</b> Lyon <b>Region</b> <b>Date</b> 08/07/1996 <b>Freight</b> 41.34
<b>Order</b> 10248 <b>Name</b> Vins et alcools Chevalier <b>City</b> Reims <b>Region</b> <b>Date</b> 04/07/1996 <b>Freight</b> 32.38		

Count : 28 Total : 1147.58

Count : 36 Total : 1571.92

Count : 16 Total : 674.93

Count : 16 Total : 674.93

Count : 16 Total : 674.93

Count : 16

8/20

Total : 674.93



# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **G**

Germany

Fri, Apr 4, 1997

<b>Order</b> 10497 <b>Name</b> Lehmanns Marktstand <b>City</b> Frankfurt a.M. <b>Region</b> <b>Date</b> 04/04/1997 <b>Freight</b> 36.21	<b>Order</b> 10488 <b>Name</b> Frankenversand <b>City</b> München <b>Region</b> <b>Date</b> 27/03/1997 <b>Freight</b> 4.93	<b>Order</b> 10468 <b>Name</b> Königlich Essen <b>City</b> Brandenburg <b>Region</b> <b>Date</b> 07/03/1997 <b>Freight</b> 44.12
<b>Order</b> 10457 <b>Name</b> Königlich Essen <b>City</b> Brandenburg <b>Region</b> <b>Date</b> 25/02/1997 <b>Freight</b> 11.57	<b>Order</b> 10456 <b>Name</b> Königlich Essen <b>City</b> Brandenburg <b>Region</b> <b>Date</b> 25/02/1997 <b>Freight</b> 8.12	<b>Order</b> 10451 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 19/02/1997 <b>Freight</b> 189.09
<b>Order</b> 10446 <b>Name</b> Toms Spezialitäten <b>City</b> Münster <b>Region</b> <b>Date</b> 14/02/1997 <b>Freight</b> 14.68	<b>Order</b> 10438 <b>Name</b> Toms Spezialitäten <b>City</b> Münster <b>Region</b> <b>Date</b> 06/02/1997 <b>Freight</b> 8.24	<b>Order</b> 10418 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 17/01/1997 <b>Freight</b> 17.55
<b>Order</b> 10407 <b>Name</b> Ottilies Käseladen <b>City</b> Köln <b>Region</b> <b>Date</b> 07/01/1997 <b>Freight</b> 91.48	<b>Order</b> 10396 <b>Name</b> Frankenversand <b>City</b> München <b>Region</b> <b>Date</b> 27/12/1996 <b>Freight</b> 135.35	<b>Order</b> 10391 <b>Name</b> Drachenblut Delikatessen <b>City</b> Aachen <b>Region</b> <b>Date</b> 23/12/1996 <b>Freight</b> 5.45
<b>Order</b> 10363 <b>Name</b> Drachenblut Delikatessen <b>City</b> Aachen <b>Region</b> <b>Date</b> 26/11/1996 <b>Freight</b> 30.54	<b>Order</b> 10361 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 22/11/1996 <b>Freight</b> 183.17	<b>Order</b> 10356 <b>Name</b> Die Wandernde Kuh <b>City</b> Stuttgart <b>Region</b> <b>Date</b> 18/11/1996 <b>Freight</b> 36.71
<b>Order</b> 10348 <b>Name</b> Die Wandernde Kuh <b>City</b> Stuttgart <b>Region</b> <b>Date</b> 07/11/1996 <b>Freight</b> 0.78	<b>Order</b> 10345 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 04/11/1996 <b>Freight</b> 249.06	<b>Order</b> 10343 <b>Name</b> Lehmanns Marktstand <b>City</b> Frankfurt a.M. <b>Region</b> <b>Date</b> 31/10/1996 <b>Freight</b> 110.37
<b>Order</b> 10342 <b>Name</b> Frankenversand <b>City</b> München <b>Region</b> <b>Date</b> 30/10/1996 <b>Freight</b> 54.83	<b>Order</b> 10337 <b>Name</b> Frankenversand <b>City</b> München <b>Region</b> <b>Date</b> 24/10/1996 <b>Freight</b> 108.26	<b>Order</b> 10325 <b>Name</b> Königlich Essen <b>City</b> Brandenburg <b>Region</b> <b>Date</b> 09/10/1996 <b>Freight</b> 64.86

Count : 21	Total : 1405.37	Count : 21	Total : 1405.37	Count : 21	Total : 1405.37
------------	-----------------	------------	-----------------	------------	-----------------

Count : 21	9/20	Total : 1405.37
------------	------	-----------------

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **G**

Germany

Fri, Apr 4, 1997

<b>Order</b> 10323 <b>Name</b> Königlich Essen <b>City</b> Brandenburg <b>Region</b> <b>Date</b> 07/10/1996 <b>Freight</b> 4.88	<b>Order</b> 10313 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 24/09/1996 <b>Freight</b> 1.96	<b>Order</b> 10312 <b>Name</b> Die Wandernde Kuh <b>City</b> Stuttgart <b>Region</b> <b>Date</b> 23/09/1996 <b>Freight</b> 40.26
<b>Order</b> 10301 <b>Name</b> Die Wandernde Kuh <b>City</b> Stuttgart <b>Region</b> <b>Date</b> 09/09/1996 <b>Freight</b> 45.08	<b>Order</b> 10286 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 21/08/1996 <b>Freight</b> 229.24	<b>Order</b> 10285 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 20/08/1996 <b>Freight</b> 76.83
<b>Order</b> 10284 <b>Name</b> Lehmanns Marktstand <b>City</b> Frankfurt a.M. <b>Region</b> <b>Date</b> 19/08/1996 <b>Freight</b> 76.56	<b>Order</b> 10279 <b>Name</b> Lehmanns Marktstand <b>City</b> Frankfurt a.M. <b>Region</b> <b>Date</b> 13/08/1996 <b>Freight</b> 25.83	<b>Order</b> 10277 <b>Name</b> Morgenstern Gesundkost <b>City</b> Leipzig <b>Region</b> <b>Date</b> 09/08/1996 <b>Freight</b> 125.77
<b>Order</b> 10273 <b>Name</b> QUICK-Stop <b>City</b> Cunewalde <b>Region</b> <b>Date</b> 05/08/1996 <b>Freight</b> 76.07	<b>Order</b> 10267 <b>Name</b> Frankenversand <b>City</b> München <b>Region</b> <b>Date</b> 29/07/1996 <b>Freight</b> 208.58	<b>Order</b> <b>10260</b> <b>Name</b> Ottilies Käseladen <b>City</b> Köln <b>Region</b> <b>Date</b> 19/07/1996 <b>Freight</b> 55.09
<b>Order</b> 10249 <b>Name</b> Toms Spezialitäten <b>City</b> Münster <b>Region</b> <b>Date</b> 05/07/1996 <b>Freight</b> 11.61		

Count : 34 Total : 2383.13

Count : 34 Total : 2383.13

Count : 13 Total : 977.76

Count : 13 Total : 977.76

Count : 13 Total : 977.76

Count : 13

10/20

Total : 977.76

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **I**

Ireland

Wed, Jan 29, 1997

<b>Order</b> 10429 <b>Name</b> Hungry Owl All-Night Grocers <b>City</b> Cork <b>Region</b> Co. Cork <b>Date</b> 29/01/1997 <b>Freight</b> 56.63	<b>Order</b> <b>10380</b> <b>Name</b> Hungry Owl All-Night Grocers <b>City</b> Cork <b>Region</b> Co. Cork <b>Date</b> 12/12/1996 <b>Freight</b> 35.03	<b>Order</b> 10373 <b>Name</b> Hungry Owl All-Night Grocers <b>City</b> Cork <b>Region</b> Co. Cork <b>Date</b> 05/12/1996 <b>Freight</b> 124.12
<b>Order</b> 10335 <b>Name</b> Hungry Owl All-Night Grocers <b>City</b> Cork <b>Region</b> Co. Cork <b>Date</b> 22/10/1996 <b>Freight</b> 42.11	<b>Order</b> 10309 <b>Name</b> Hungry Owl All-Night Grocers <b>City</b> Cork <b>Region</b> Co. Cork <b>Date</b> 19/09/1996 <b>Freight</b> 47.30	<b>Order</b> 10298 <b>Name</b> Hungry Owl All-Night Grocers <b>City</b> Cork <b>Region</b> Co. Cork <b>Date</b> 05/09/1996 <b>Freight</b> 168.22

Count : 6 Total : 473.41

Italy

Thu, Mar 6, 1997

<b>Order</b> 10467 <b>Name</b> Magazzini Alimentari Riuniti <b>City</b> Bergamo <b>Region</b> <b>Date</b> 06/03/1997 <b>Freight</b> 4.93	<b>Order</b> 10443 <b>Name</b> Reggiani Caseifici <b>City</b> Reggio Emilia <b>Region</b> <b>Date</b> 12/02/1997 <b>Freight</b> 13.95	<b>Order</b> 10428 <b>Name</b> Reggiani Caseifici <b>City</b> Reggio Emilia <b>Region</b> <b>Date</b> 28/01/1997 <b>Freight</b> 11.09
<b>Order</b> 10422 <b>Name</b> Franchi S.p.A. <b>City</b> Torino <b>Region</b> <b>Date</b> 22/01/1997 <b>Freight</b> 3.02	<b>Order</b> 10404 <b>Name</b> Magazzini Alimentari Riuniti <b>City</b> Bergamo <b>Region</b> <b>Date</b> 03/01/1997 <b>Freight</b> 155.97	<b>Order</b> <b>10300</b> <b>Name</b> Magazzini Alimentari Riuniti <b>City</b> Bergamo <b>Region</b> <b>Date</b> 09/09/1996 <b>Freight</b> 17.68
<b>Order</b> 10288 <b>Name</b> Reggiani Caseifici <b>City</b> Reggio Emilia <b>Region</b> <b>Date</b> 23/08/1996 <b>Freight</b> 7.45	<b>Order</b> 10275 <b>Name</b> Magazzini Alimentari Riuniti <b>City</b> Bergamo <b>Region</b> <b>Date</b> 07/08/1996 <b>Freight</b> 26.93	

Count : 8 Total : 241.02

Count : 14 Total : 714.43

Count : 14 Total : 714.43

Count : 14 Total : 714.43

Count : 14 Total : 714.43

Count : 14 11/20 Total : 714.43

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **M**

**Mexico**

**Thu, Mar 13, 1997**

<b>Order</b> 10474 <b>Name</b> Pericles Comidas clásicas <b>City</b> México D.F. <b>Region</b> <b>Date</b> 13/03/1997 <b>Freight</b> 83.49	<b>Order</b> 10365 <b>Name</b> Antonio Moreno Taquería <b>City</b> México D.F. <b>Region</b> <b>Date</b> 27/11/1996 <b>Freight</b> 22.00	<b>Order</b> 10354 <b>Name</b> Pericles Comidas clásicas <b>City</b> México D.F. <b>Region</b> <b>Date</b> 14/11/1996 <b>Freight</b> 53.80
<b>Order</b> 10322 <b>Name</b> Pericles Comidas clásicas <b>City</b> México D.F. <b>Region</b> <b>Date</b> 04/10/1996 <b>Freight</b> 0.40	<b>Order</b> 10319 <b>Name</b> Tortuga Restaurante <b>City</b> México D.F. <b>Region</b> <b>Date</b> 02/10/1996 <b>Freight</b> 64.50	<b>Order</b> 10308 <b>Name</b> Ana Trujillo Emparedados y helados <b>City</b> México D.F. <b>Region</b> <b>Date</b> 18/09/1996 <b>Freight</b> 1.61
<b>Order</b> 10304 <b>Name</b> Tortuga Restaurante <b>City</b> México D.F. <b>Region</b> <b>Date</b> 12/09/1996 <b>Freight</b> 63.79	<b>Order</b> 10293 <b>Name</b> Tortuga Restaurante <b>City</b> México D.F. <b>Region</b> <b>Date</b> 29/08/1996 <b>Freight</b> 21.18	<b>Order</b> 10276 <b>Name</b> Tortuga Restaurante <b>City</b> México D.F. <b>Region</b> <b>Date</b> 08/08/1996 <b>Freight</b> 13.84
<b>Order</b> 10259 <b>Name</b> Centro comercial Moctezuma <b>City</b> México D.F. <b>Region</b> <b>Date</b> 18/07/1996 <b>Freight</b> 3.25		

**Count : 10      Total : 327.86**

**Count : 10      Total : 327.86**

**Count : 10      Total : 327.86      Count : 10      Total : 327.86      Count : 10      Total : 327.86**

**Count : 10      12/20      Total : 327.86**

Northwind Order List

Max order ID is : 10500

OrdersOrdersOrders

Countries Starting With Letter : N

NorwayWed, Dec 18, 1996

Order	10387
Name	Santé Gourmet
City	Stavern
Region	
Date	18/12/1996
Freight	93.63

Count : 1Total : 93.63

Count : 1Total : 93.63

Count : 1Total : 93.63Count : 1Total : 93.63Count : 1Total : 93.63

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **P**

**Poland**

**Thu, Dec 5, 1996**

<b>Order</b>	10374
<b>Name</b>	Wolski Zajazd
<b>City</b>	Warszawa
<b>Region</b>	
<b>Date</b>	05/12/1996
<b>Freight</b>	3.94

**Count :** 1 **Total :** 3.94

**Portugal**

**Mon, Mar 31, 1997**

<b>Order</b>	10491
<b>Name</b>	Furia Bacalhau e Frutos do Mar
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	31/03/1997
<b>Freight</b>	16.96

<b>Order</b>	10477
<b>Name</b>	Princesa Isabel Vinhos
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	17/03/1997
<b>Freight</b>	13.02

<b>Order</b>	10464
<b>Name</b>	Furia Bacalhau e Frutos do Mar
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	04/03/1997
<b>Freight</b>	89.00

<b>Order</b>	10433
<b>Name</b>	Princesa Isabel Vinhos
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	03/02/1997
<b>Freight</b>	73.83

<b>Order</b>	10397
<b>Name</b>	Princesa Isabel Vinhos
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	27/12/1996
<b>Freight</b>	60.26

<b>Order</b>	10352
<b>Name</b>	Furia Bacalhau e Frutos do Mar
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	12/11/1996
<b>Freight</b>	1.30

<b>Order</b>	10336
<b>Name</b>	Princesa Isabel Vinhos
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	23/10/1996
<b>Freight</b>	15.51

<b>Order</b>	10328
<b>Name</b>	Furia Bacalhau e Frutos do Mar
<b>City</b>	Lisboa
<b>Region</b>	
<b>Date</b>	14/10/1996
<b>Freight</b>	87.03

**Count :** 8 **Total :** 356.91

**Count :** 9 **Total :** 360.85

**Count :** 9 **Total :** 360.85

**Count :** 9 **Total :** 360.85

**Count :** 9 **Total :** 360.85

**Count :** 9 **14/20** **Total :** 360.85

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **S**

Spain

Mon, Jan 27, 1997

<b>Order</b> 10426 <b>Name</b> Galería del gastrónomo <b>City</b> Barcelona <b>Region</b> <b>Date</b> 27/01/1997 <b>Freight</b> 18.69	<b>Order</b> 10366 <b>Name</b> Galería del gastrónomo <b>City</b> Barcelona <b>Region</b> <b>Date</b> 28/11/1996 <b>Freight</b> 10.14	<b>Order</b> 10326 <b>Name</b> Bólido Comidas preparadas <b>City</b> Madrid <b>Region</b> <b>Date</b> 10/10/1996 <b>Freight</b> 77.92
<b>Order</b> 10306 <b>Name</b> Romero y tomillo <b>City</b> Madrid <b>Region</b> <b>Date</b> 16/09/1996 <b>Freight</b> 7.56	<b>Order</b> 10303 <b>Name</b> Godos Cocina Típica <b>City</b> Sevilla <b>Region</b> <b>Date</b> 11/09/1996 <b>Freight</b> 107.83	<b>Order</b> 10282 <b>Name</b> Romero y tomillo <b>City</b> Madrid <b>Region</b> <b>Date</b> 15/08/1996 <b>Freight</b> 12.69
<b>Order</b> 10281 <b>Name</b> Romero y tomillo <b>City</b> Madrid <b>Region</b> <b>Date</b> 14/08/1996 <b>Freight</b> 2.94		
<b>Count :</b> 7 <b>Total :</b> 237.77		

Sweden

Fri, Feb 28, 1997

<b>Order</b> 10460 <b>Name</b> Folk och få HB <b>City</b> Bräcke <b>Region</b> <b>Date</b> 28/02/1997 <b>Freight</b> 16.27	<b>Order</b> 10445 <b>Name</b> Berglunds snabbköp <b>City</b> Luleå <b>Region</b> <b>Date</b> 13/02/1997 <b>Freight</b> 9.30	<b>Order</b> 10444 <b>Name</b> Berglunds snabbköp <b>City</b> Luleå <b>Region</b> <b>Date</b> 12/02/1997 <b>Freight</b> 3.50
<b>Order</b> 10434 <b>Name</b> Folk och få HB <b>City</b> Bräcke <b>Region</b> <b>Date</b> 03/02/1997 <b>Freight</b> 17.92	<b>Order</b> 10384 <b>Name</b> Berglunds snabbköp <b>City</b> Luleå <b>Region</b> <b>Date</b> 16/12/1996 <b>Freight</b> 168.64	<b>Order</b> 10378 <b>Name</b> Folk och få HB <b>City</b> Bräcke <b>Region</b> <b>Date</b> 10/12/1996 <b>Freight</b> 5.44
<b>Order</b> 10327 <b>Name</b> Folk och få HB <b>City</b> Bräcke <b>Region</b> <b>Date</b> 11/10/1996 <b>Freight</b> 63.36	<b>Order</b> 10280 <b>Name</b> Berglunds snabbköp <b>City</b> Luleå <b>Region</b> <b>Date</b> 14/08/1996 <b>Freight</b> 8.98	<b>Order</b> 10278 <b>Name</b> Berglunds snabbköp <b>City</b> Luleå <b>Region</b> <b>Date</b> 12/08/1996 <b>Freight</b> 92.69
<b>Order</b> 10264 <b>Name</b> Folk och få HB <b>City</b> Bräcke <b>Region</b> <b>Date</b> 24/07/1996 <b>Freight</b> 3.67		
<b>Count :</b> 10 <b>Total :</b> 389.77		

Count : 17 Total : 627.54

Count : 17 Total : 627.54

Count : 17 Total : 627.54

Count : 17 15/20 Total : 627.54

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **S**

Switzerland

Mon, Jan 20, 1997

Order	10419
Name	Richter Supermarkt
City	Genève
Region	
Date	20/01/1997
Freight	137.35

Order	<b>10370</b>
Name	Chop-suey Chinese
City	Bern
Region	
Date	03/12/1996
Freight	1.17

Order	10255
Name	Richter Supermarkt
City	Genève
Region	
Date	12/07/1996
Freight	148.33

Order	10254
Name	Chop-suey Chinese
City	Bern
Region	
Date	11/07/1996
Freight	22.98

Count :	4	Total :	309.83
---------	---	---------	--------

Count :	21	Total :	937.37
---------	----	---------	--------

Count :	4	Total :	309.83	Count :	4	Total :	309.83	Count :	4	Total :	309.83
---------	---	---------	--------	---------	---	---------	--------	---------	---	---------	--------

Count :	4	16/20	Total :	309.83
---------	---	-------	---------	--------



# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **U**

**UK** **Mon, Mar 24, 1997**

<b>Order</b> 10484 <b>Name</b> B's Beverages <b>City</b> London <b>Region</b> <b>Date</b> 24/03/1997 <b>Freight</b> 6.88	<b>Order</b> 10473 <b>Name</b> Island Trading <b>City</b> Cowes <b>Region</b> Isle of Wight <b>Date</b> 13/03/1997 <b>Freight</b> 16.37	<b>Order</b> 10472 <b>Name</b> Seven Seas Imports <b>City</b> London <b>Region</b> <b>Date</b> 12/03/1997 <b>Freight</b> 4.20
<b>Order</b> 10471 <b>Name</b> B's Beverages <b>City</b> London <b>Region</b> <b>Date</b> 11/03/1997 <b>Freight</b> 45.59	<b>Order</b> 10462 <b>Name</b> Consolidated Holdings <b>City</b> London <b>Region</b> <b>Date</b> 03/03/1997 <b>Freight</b> 6.17	<b>Order</b> 10453 <b>Name</b> Around the Horn <b>City</b> Colchester <b>Region</b> Essex <b>Date</b> 21/02/1997 <b>Freight</b> 25.36
<b>Order</b> 10435 <b>Name</b> Consolidated Holdings <b>City</b> London <b>Region</b> <b>Date</b> 04/02/1997 <b>Freight</b> 9.21	<b>Order</b> <b>10400</b> <b>Name</b> Eastern Connection <b>City</b> London <b>Region</b> <b>Date</b> 01/01/1997 <b>Freight</b> 83.93	<b>Order</b> 10388 <b>Name</b> Seven Seas Imports <b>City</b> London <b>Region</b> <b>Date</b> 19/12/1996 <b>Freight</b> 34.86
<b>Order</b> 10383 <b>Name</b> Around the Horn <b>City</b> Colchester <b>Region</b> Essex <b>Date</b> 16/12/1996 <b>Freight</b> 34.24	<b>Order</b> 10377 <b>Name</b> Seven Seas Imports <b>City</b> London <b>Region</b> <b>Date</b> 09/12/1996 <b>Freight</b> 22.21	<b>Order</b> 10364 <b>Name</b> Eastern Connection <b>City</b> London <b>Region</b> <b>Date</b> 26/11/1996 <b>Freight</b> 71.97
<b>Order</b> 10359 <b>Name</b> Seven Seas Imports <b>City</b> London <b>Region</b> <b>Date</b> 21/11/1996 <b>Freight</b> 288.43	<b>Order</b> 10355 <b>Name</b> Around the Horn <b>City</b> Colchester <b>Region</b> Essex <b>Date</b> 15/11/1996 <b>Freight</b> 41.95	<b>Order</b> 10321 <b>Name</b> Island Trading <b>City</b> Cowes <b>Region</b> Isle of Wight <b>Date</b> 03/10/1996 <b>Freight</b> 3.43
<b>Order</b> 10318 <b>Name</b> Island Trading <b>City</b> Cowes <b>Region</b> Isle of Wight <b>Date</b> 01/10/1996 <b>Freight</b> 4.73	<b>Order</b> 10315 <b>Name</b> Island Trading <b>City</b> Cowes <b>Region</b> Isle of Wight <b>Date</b> 26/09/1996 <b>Freight</b> 41.76	<b>Order</b> 10289 <b>Name</b> B's Beverages <b>City</b> London <b>Region</b> <b>Date</b> 26/08/1996 <b>Freight</b> 22.77

**Count : 18 Total : 764.06**

**USA** **Mon, Mar 24, 1997**

<b>Order</b> 10483 <b>Name</b> White Clover Markets <b>City</b> Seattle <b>Region</b> WA <b>Date</b> 24/03/1997 <b>Freight</b> 15.28	<b>Order</b> 10482 <b>Name</b> Lazy K Kountry Store <b>City</b> Walla Walla <b>Region</b> WA <b>Date</b> 21/03/1997 <b>Freight</b> 7.48	<b>Order</b> 10479 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 19/03/1997 <b>Freight</b> 708.95
---	--	--

**Count : 21 Total : 1495.77** **Count : 21 Total : 1495.77** **Count : 21 Total : 1495.77**

**Count : 21 17/20 Total : 1495.77**

# Northwind Order List

Max order ID is : 10500

Orders Orders Orders

Countries Starting With Letter : **U**

USA

Mon, Mar 24, 1997

<b>Order</b> 10469 <b>Name</b> White Clover Markets <b>City</b> Seattle <b>Region</b> WA <b>Date</b> 10/03/1997 <b>Freight</b> 60.18	<b>Order</b> 10452 <b>Name</b> Save-a-lot Markets <b>City</b> Boise <b>Region</b> ID <b>Date</b> 20/02/1997 <b>Freight</b> 140.26	<b>Order</b> 10441 <b>Name</b> Old World Delicatessen <b>City</b> Anchorage <b>Region</b> AK <b>Date</b> 10/02/1997 <b>Freight</b> 73.02
<b>Order</b> <b>10440</b> <b>Name</b> Save-a-lot Markets <b>City</b> Boise <b>Region</b> ID <b>Date</b> 10/02/1997 <b>Freight</b> 86.53	<b>Order</b> 10432 <b>Name</b> Split Rail Beer & Ale <b>City</b> Lander <b>Region</b> WY <b>Date</b> 31/01/1997 <b>Freight</b> 4.34	<b>Order</b> 10415 <b>Name</b> Hungry Coyote Import Store <b>City</b> Elgin <b>Region</b> OR <b>Date</b> 15/01/1997 <b>Freight</b> 0.20
<b>Order</b> 10401 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 01/01/1997 <b>Freight</b> 12.51	<b>Order</b> 10398 <b>Name</b> Save-a-lot Markets <b>City</b> Boise <b>Region</b> ID <b>Date</b> 30/12/1996 <b>Freight</b> 89.16	<b>Order</b> 10394 <b>Name</b> Hungry Coyote Import Store <b>City</b> Elgin <b>Region</b> OR <b>Date</b> 25/12/1996 <b>Freight</b> 30.34
<b>Order</b> 10393 <b>Name</b> Save-a-lot Markets <b>City</b> Boise <b>Region</b> ID <b>Date</b> 25/12/1996 <b>Freight</b> 126.56	<b>Order</b> 10385 <b>Name</b> Split Rail Beer & Ale <b>City</b> Lander <b>Region</b> WY <b>Date</b> 17/12/1996 <b>Freight</b> 30.96	<b>Order</b> 10375 <b>Name</b> Hungry Coyote Import Store <b>City</b> Elgin <b>Region</b> OR <b>Date</b> 06/12/1996 <b>Freight</b> 20.12
<b>Order</b> 10369 <b>Name</b> Split Rail Beer & Ale <b>City</b> Lander <b>Region</b> WY <b>Date</b> 02/12/1996 <b>Freight</b> 195.68	<b>Order</b> 10349 <b>Name</b> Split Rail Beer & Ale <b>City</b> Lander <b>Region</b> WY <b>Date</b> 08/11/1996 <b>Freight</b> 8.63	<b>Order</b> 10346 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 05/11/1996 <b>Freight</b> 142.08
<b>Order</b> 10344 <b>Name</b> White Clover Markets <b>City</b> Seattle <b>Region</b> WA <b>Date</b> 01/11/1996 <b>Freight</b> 23.29	<b>Order</b> 10338 <b>Name</b> Old World Delicatessen <b>City</b> Anchorage <b>Region</b> AK <b>Date</b> 25/10/1996 <b>Freight</b> 84.21	<b>Order</b> 10329 <b>Name</b> Split Rail Beer & Ale <b>City</b> Lander <b>Region</b> WY <b>Date</b> 15/10/1996 <b>Freight</b> 191.67
<b>Order</b> 10324 <b>Name</b> Save-a-lot Markets <b>City</b> Boise <b>Region</b> ID <b>Date</b> 08/10/1996 <b>Freight</b> 214.27	<b>Order</b> 10317 <b>Name</b> Lonesome Pine Restaurant <b>City</b> Portland <b>Region</b> OR <b>Date</b> 30/09/1996 <b>Freight</b> 12.69	<b>Order</b> 10316 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 27/09/1996 <b>Freight</b> 150.15

Count : 21 Total : 1696.85 Count : 21 Total : 1696.85 Count : 21 Total : 1696.85

Count : 21 18/20 Total : 1696.85

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **U**

**USA**

**Mon, Mar 24, 1997**

<b>Order</b> 10314 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 25/09/1996 <b>Freight</b> 74.16	<b>Order</b> <b>10310</b> <b>Name</b> The Big Cheese <b>City</b> Portland <b>Region</b> OR <b>Date</b> 20/09/1996 <b>Freight</b> 17.52	<b>Order</b> 10307 <b>Name</b> Lonesome Pine Restaurant <b>City</b> Portland <b>Region</b> OR <b>Date</b> 17/09/1996 <b>Freight</b> 0.56
<b>Order</b> 10305 <b>Name</b> Old World Delicatessen <b>City</b> Anchorage <b>Region</b> AK <b>Date</b> 13/09/1996 <b>Freight</b> 257.62	<b>Order</b> 10294 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 30/08/1996 <b>Freight</b> 147.26	<b>Order</b> 10272 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 02/08/1996 <b>Freight</b> 98.03
<b>Order</b> 10271 <b>Name</b> Split Rail Beer & Ale <b>City</b> Lander <b>Region</b> WY <b>Date</b> 01/08/1996 <b>Freight</b> 4.54	<b>Order</b> 10269 <b>Name</b> White Clover Markets <b>City</b> Seattle <b>Region</b> WA <b>Date</b> 31/07/1996 <b>Freight</b> 4.56	<b>Order</b> 10262 <b>Name</b> Rattlesnake Canyon Grocery <b>City</b> Albuquerque <b>Region</b> NM <b>Date</b> 22/07/1996 <b>Freight</b> 48.29

**Count : 33      Total : 3081.10**

**Count : 51      Total : 3845.16**

**Count : 9      Total : 652.54      Count : 9      Total : 652.54      Count : 9      Total : 652.54**

**Count : 9      19/20      Total : 652.54**

# Northwind Order List

Max order ID is : 10500

Orders	Orders	Orders
--------	--------	--------

Countries Starting With Letter : **V**

Venezuela

Tue, Apr 8, 1997

<b>Order</b> 10499 <b>Name</b> LILA-Supermercado <b>City</b> Barquisimeto <b>Region</b> Lara <b>Date</b> 08/04/1997 <b>Freight</b> 102.02	<b>Order</b> 10498 <b>Name</b> HILARION-Abastos <b>City</b> San Cristóbal <b>Region</b> Táchira <b>Date</b> 07/04/1997 <b>Freight</b> 29.75	<b>Order</b> <b>10490</b> <b>Name</b> HILARION-Abastos <b>City</b> San Cristóbal <b>Region</b> Táchira <b>Date</b> 31/03/1997 <b>Freight</b> 210.19
<b>Order</b> 10486 <b>Name</b> HILARION-Abastos <b>City</b> San Cristóbal <b>Region</b> Táchira <b>Date</b> 26/03/1997 <b>Freight</b> 30.53	<b>Order</b> 10485 <b>Name</b> LINO-Delicateses <b>City</b> I. de Margarita <b>Region</b> Nueva Esparta <b>Date</b> 25/03/1997 <b>Freight</b> 64.45	<b>Order</b> 10476 <b>Name</b> HILARION-Abastos <b>City</b> San Cristóbal <b>Region</b> Táchira <b>Date</b> 17/03/1997 <b>Freight</b> 4.41
<b>Order</b> 10461 <b>Name</b> LILA-Supermercado <b>City</b> Barquisimeto <b>Region</b> Lara <b>Date</b> 28/02/1997 <b>Freight</b> 148.61	<b>Order</b> 10405 <b>Name</b> LINO-Delicateses <b>City</b> I. de Margarita <b>Region</b> Nueva Esparta <b>Date</b> 06/01/1997 <b>Freight</b> 34.82	<b>Order</b> 10395 <b>Name</b> HILARION-Abastos <b>City</b> San Cristóbal <b>Region</b> Táchira <b>Date</b> 26/12/1996 <b>Freight</b> 184.41
<b>Order</b> 10381 <b>Name</b> LILA-Supermercado <b>City</b> Barquisimeto <b>Region</b> Lara <b>Date</b> 12/12/1996 <b>Freight</b> 7.99	<b>Order</b> 10357 <b>Name</b> LILA-Supermercado <b>City</b> Barquisimeto <b>Region</b> Lara <b>Date</b> 19/11/1996 <b>Freight</b> 34.88	<b>Order</b> <b>10330</b> <b>Name</b> LILA-Supermercado <b>City</b> Barquisimeto <b>Region</b> Lara <b>Date</b> 16/10/1996 <b>Freight</b> 12.75
<b>Order</b> 10296 <b>Name</b> LILA-Supermercado <b>City</b> Barquisimeto <b>Region</b> Lara <b>Date</b> 03/09/1996 <b>Freight</b> 0.12	<b>Order</b> 10283 <b>Name</b> LILA-Supermercado <b>City</b> Barquisimeto <b>Region</b> Lara <b>Date</b> 16/08/1996 <b>Freight</b> 84.81	<b>Order</b> 10268 <b>Name</b> GROSELLA-Restaurante <b>City</b> Caracas <b>Region</b> DF <b>Date</b> 30/07/1996 <b>Freight</b> 66.29
<b>Order</b> 10257 <b>Name</b> HILARION-Abastos <b>City</b> San Cristóbal <b>Region</b> Táchira <b>Date</b> 16/07/1996 <b>Freight</b> 81.91		

Count : 16 Total : 1097.94

Count : 16 Total : 1097.94



That's All Folks!  
Hei\_remind\_me\_to\_put myself up for  
abduction. END!

Count : 16 Total : 1097.94

Count : 16 Total : 1097.94

Count : 16 Total : 1097.94

Count : 16

20/20

Total : 1097.94