

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

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
-------	------------	------	---------

Countries Starting With Letter : **A**

Argentina		Mon, Feb 17, 1997	
10448	Rancho grande, Buenos Aires	17/02/1997	38.82
10409	Océano Atlántico Ltda., Buenos Aires	09/01/1997	29.83
Count : 2		Total :	68.65
Austria		Fri, Mar 28, 1997	
10489	Piccolo und mehr, Salzburg	28/03/1997	5.29
10442	Ernst Handel, Graz	11/02/1997	47.94
10430	Ernst Handel, Graz	30/01/1997	458.78
10427	Piccolo und mehr, Salzburg	27/01/1997	31.29
10403	Ernst Handel, Graz	03/01/1997	73.79
10402	Ernst Handel, Graz	02/01/1997	67.88
10392	Piccolo und mehr, Salzburg	24/12/1996	122.46
10390	Ernst Handel, Graz	23/12/1996	126.38
10382	Ernst Handel, Graz	13/12/1996	94.77
10368	Ernst Handel, Graz	29/11/1996	101.95
10353	Piccolo und mehr, Salzburg	13/11/1996	360.63
10351	Ernst Handel, Graz	11/11/1996	162.33
10263	Ernst Handel, Graz	23/07/1996	146.06
10258	Ernst Handel, Graz	17/07/1996	140.51
Count : 14		Total :	1940.06
Count : 16		Total :	2008.71

Order	Name, City	Date	Freight
-------	------------	------	---------

Countries Starting With Letter : **B**

Belgium		Fri, Mar 14, 1997	
10475	Suprêmes délices, Charleroi	14/03/1997	68.52
10463	Suprêmes délices, Charleroi	04/03/1997	14.78
10458	Suprêmes délices, Charleroi	26/02/1997	147.06
10302	Suprêmes délices, Charleroi	10/09/1996	6.27
10252	Suprêmes délices, Charleroi	09/07/1996	51.30
Count : 5		Total :	287.93
Brazil		Fri, Apr 4, 1997	
10496	Tradição Hipermercados, Sao Paulo	04/04/1997	46.77
10494	Comércio Mineiro, Sao Paulo	02/04/1997	65.99
10487	Queen Cozinha, Sao Paulo	26/03/1997	71.07
10481	Ricardo Adocicados, Rio de Janeiro	20/03/1997	64.33
10466	Comércio Mineiro, Sao Paulo	06/03/1997	11.93
10447	Ricardo Adocicados, Rio de Janeiro	14/02/1997	68.66
10423	Gourmet Lanchonetes, Campinas	23/01/1997	24.50
10421	Que Delícia, Rio de Janeiro	21/01/1997	99.23
10420	Wellington Importadora, Resende	21/01/1997	44.12
10414	Familia Arquibaldo, Sao Paulo	14/01/1997	21.48
10406	Queen Cozinha, Sao Paulo	07/01/1997	108.04
10386	Familia Arquibaldo, Sao Paulo	18/12/1996	13.99
10379	Que Delícia, Rio de Janeiro	11/12/1996	45.03
10372	Queen Cozinha, Sao Paulo	04/12/1996	890.78
10347	Familia Arquibaldo, Sao Paulo	06/11/1996	3.10
10299	Ricardo Adocicados, Rio de Janeiro	06/09/1996	29.76
10292	Tradição Hipermercados, Sao Paulo	28/08/1996	1.35
10291	Que Delícia, Rio de Janeiro	27/08/1996	6.40
10290	Comércio Mineiro, Sao Paulo	27/08/1996	79.70
10287	Ricardo Adocicados, Rio de Janeiro	22/08/1996	12.76
10261	Que Delícia, Rio de Janeiro	19/07/1996	3.05
10256	Wellington Importadora, Resende	15/07/1996	13.97
10253	Hanari Carnes, Rio de Janeiro	10/07/1996	58.17
10250	Hanari Carnes, Rio de Janeiro	08/07/1996	65.83
Count : 29		Total :	2137.94

Count : 16	Total : 2008.71
-------------------	------------------------

Count : 45	1/8	Total : 4146.65
-------------------	------------	------------------------

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
-------	------------	------	---------

Countries Starting With Letter : **B**

Brazil		Fri, Apr 4, 1997	
Count :	24	Total :	1850.01
Count :	29	Total :	2137.94

Order	Name, City	Date	Freight
-------	------------	------	---------

Countries Starting With Letter : **C**

Canada		Thu, Apr 3, 1997		
10495	Laughing Bacchus Wine Cellars, Vancouver	BC	03/04/1997	4.65
10492	Bottom-Dollar Markets, Tsawassen	BC	01/04/1997	62.89
10439	Mère Paillarde, Montréal	Québe	07/02/1997	4.07
10431	Bottom-Dollar Markets, Tsawassen	BC	30/01/1997	44.17
10424	Mère Paillarde, Montréal	Québe	23/01/1997	370.61
10411	Bottom-Dollar Markets, Tsawassen	BC	10/01/1997	23.65
10410	Bottom-Dollar Markets, Tsawassen	BC	10/01/1997	2.40
10389	Bottom-Dollar Markets, Tsawassen	BC	20/12/1996	47.42
10376	Mère Paillarde, Montréal	Québe	09/12/1996	20.39
10339	Mère Paillarde, Montréal	Québe	28/10/1996	15.66
10332	Mère Paillarde, Montréal	Québe	17/10/1996	52.84
Count : 11		Total :		648.75
Count : 11		Total :		648.75

Count :	29	Total :	2137.94
----------------	-----------	----------------	----------------

Count :	11	Total :	648.75
----------------	-----------	----------------	---------------

Count :	11	2/8	Total :	648.75
----------------	-----------	------------	----------------	---------------

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
Countries Starting With Letter : D			
Denmark		Wed, Mar 5, 1997	
10465	Vaffeljernet, Århus	05/03/1997	145.04
10417	Simons bistro, Kobenhavn	16/01/1997	70.29
10399	Vaffeljernet, Århus	31/12/1996	27.36
10367	Vaffeljernet, Århus	28/11/1996	13.55
10341	Simons bistro, Kobenhavn	29/10/1996	26.78
Count :	5	Total :	283.02
Count :	5	Total :	283.02

Order	Name, City	Date	Freight
Countries Starting With Letter : F			
Finland		Mon, Feb 24, 1997	
10455	Wartian Herkku, Oulu	24/02/1997	180.45
10437	Wartian Herkku, Oulu	05/02/1997	19.97
10416	Wartian Herkku, Oulu	16/01/1997	22.72
10412	Wartian Herkku, Oulu	13/01/1997	3.77
10333	Wartian Herkku, Oulu	18/10/1996	0.59
10320	Wartian Herkku, Oulu	03/10/1996	34.57
10270	Wartian Herkku, Oulu	01/08/1996	136.54

Count :	8	Total :	424.34
---------	---	---------	--------

France Wed, Apr 9, 1997

10500	La maison d'Asie, Toulouse	09/04/1997	42.68
10493	La maison d'Asie, Toulouse	02/04/1997	10.64
10480	Folies gourmandes, Lille	20/03/1997	1.35
10478	Victuailles en stock, Lyon	18/03/1997	4.81
10470	Bon app', Marseille	11/03/1997	64.56

10459	Victuailles en stock, Lyon	27/02/1997	25.09
10454	La maison d'Asie, Toulouse	21/02/1997	2.74
10450	Victuailles en stock, Lyon	19/02/1997	7.23
10449	Blondel père et fils, Strasbourg	18/02/1997	53.30
10436	Blondel père et fils, Strasbourg	05/02/1997	156.66

10425	La maison d'Asie, Toulouse	24/01/1997	7.93
10413	La maison d'Asie, Toulouse	14/01/1997	95.66
10408	Folies gourmandes, Lille	08/01/1997	11.26
10371	La maison d'Asie, Toulouse	03/12/1996	0.45
10362	Bon app', Marseille	25/11/1996	96.04

10360	Blondel père et fils, Strasbourg	22/11/1996	131.70
10358	La maison d'Asie, Toulouse	20/11/1996	19.64
10350	La maison d'Asie, Toulouse	11/11/1996	64.19
10340	Bon app', Marseille	29/10/1996	166.31
10334	Victuailles en stock, Lyon	21/10/1996	8.56

10331	Bon app', Marseille	16/10/1996	10.19
10311	Du monde entier, Nantes	20/09/1996	24.69
10297	Blondel père et fils, Strasbourg	04/09/1996	5.74
10295	Vins et alcools Chevalier, Reims	02/09/1996	1.15
10274	Vins et alcools Chevalier, Reims	06/08/1996	6.01

10265	Blondel père et fils, Strasbourg	25/07/1996	55.28
10251	Victuailles en stock, Lyon	08/07/1996	41.34
10248	Vins et alcools Chevalier, Reims	04/07/1996	32.38

Count :	28	Total :	1147.58
---------	----	---------	---------

Count :	36	Total :	1571.92
---------	----	---------	---------

Count :	5	Total :	283.02
---------	---	---------	--------

Count :	36	Total :	1571.92
---------	----	---------	---------

Count :	41	3/8	Total :	1854.94
---------	----	-----	---------	---------

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
Countries Starting With Letter : G			
Germany		Fri, Apr 4, 1997	
10497	Lehmanns Marktstand, Frankfurt a.M.	04/04/1997	36.21
10488	Frankenversand, München	27/03/1997	4.93
10468	Königlich Essen, Brandenburg	07/03/1997	44.12
10457	Königlich Essen, Brandenburg	25/02/1997	11.57
10456	Königlich Essen, Brandenburg	25/02/1997	8.12
10451	QUICK-Stop, Cunewalde	19/02/1997	189.09
10446	Toms Spezialitäten, Münster	14/02/1997	14.68
10438	Toms Spezialitäten, Münster	06/02/1997	8.24
10418	QUICK-Stop, Cunewalde	17/01/1997	17.55
10407	Ottilies Käseladen, Köln	07/01/1997	91.48
10396	Frankenversand, München	27/12/1996	135.35
10391	Drachenblut Delikatessen, Aachen	23/12/1996	5.45
10363	Drachenblut Delikatessen, Aachen	26/11/1996	30.54
10361	QUICK-Stop, Cunewalde	22/11/1996	183.17
10356	Die Wandernde Kuh, Stuttgart	18/11/1996	36.71
10348	Die Wandernde Kuh, Stuttgart	07/11/1996	0.78
10345	QUICK-Stop, Cunewalde	04/11/1996	249.06
10343	Lehmanns Marktstand, Frankfurt a.M.	31/10/1996	110.37
10342	Frankenversand, München	30/10/1996	54.83
10337	Frankenversand, München	24/10/1996	108.26
10325	Königlich Essen, Brandenburg	09/10/1996	64.86
10323	Königlich Essen, Brandenburg	07/10/1996	4.88
10313	QUICK-Stop, Cunewalde	24/09/1996	1.96
10312	Die Wandernde Kuh, Stuttgart	23/09/1996	40.26
10301	Die Wandernde Kuh, Stuttgart	09/09/1996	45.08
10286	QUICK-Stop, Cunewalde	21/08/1996	229.24
10285	QUICK-Stop, Cunewalde	20/08/1996	76.83
10284	Lehmanns Marktstand, Frankfurt a.M.	19/08/1996	76.56
10279	Lehmanns Marktstand, Frankfurt a.M.	13/08/1996	25.83
10277	Morgenstern Gesundkost, Leipzig	09/08/1996	125.77
10273	QUICK-Stop, Cunewalde	05/08/1996	76.07
10267	Frankenversand, München	29/07/1996	208.58
10260	Ottilies Käseladen, Köln	19/07/1996	55.09
10249	Toms Spezialitäten, Münster	05/07/1996	11.61
Count :	34	Total :	2383.13
Count :	34	Total :	2383.13

Order	Name, City	Date	Freight
Countries Starting With Letter : I			
Ireland		Wed, Jan 29, 1997	
10429	Hungry Owl All-Night Grocers, Co. Cork	29/01/1997	56.63
10380	Hungry Owl All-Night Grocers, Co. Cork	12/12/1996	35.03
10373	Hungry Owl All-Night Grocers, Co. Cork	05/12/1996	124.12
10335	Hungry Owl All-Night Grocers, Co. Cork	22/10/1996	42.11
10309	Hungry Owl All-Night Grocers, Co. Cork	19/09/1996	47.30
10298	Hungry Owl All-Night Grocers, Co. Cork	05/09/1996	168.22
Count :	6	Total :	473.41
Italy		Thu, Mar 6, 1997	
10467	Magazzini Alimentari Riuniti, Bergamo	06/03/1997	4.93
10443	Reggiani Caseifici, Reggio Emilia	12/02/1997	13.95
10428	Reggiani Caseifici, Reggio Emilia	28/01/1997	11.09
10422	Franchi S.p.A., Torino	22/01/1997	3.02
10404	Magazzini Alimentari Riuniti, Bergamo	03/01/1997	155.97
10300	Magazzini Alimentari Riuniti, Bergamo	09/09/1996	17.68
10288	Reggiani Caseifici, Reggio Emilia	23/08/1996	7.45
10275	Magazzini Alimentari Riuniti, Bergamo	07/08/1996	26.93
Count :	8	Total :	241.02
Count :	14	Total :	714.43

Count :	34	Total :	2383.13
----------------	-----------	----------------	----------------

Count :	14	Total :	714.43
----------------	-----------	----------------	---------------

Count : 48 4/8 Total : 3097.56

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
<i>Countries Starting With Letter : M</i>			
Mexico		Thu, Mar 13, 1997	
10474	Pericles Comidas clásicas, México D.F.	13/03/1997	83.49
10365	Antonio Moreno Taquería, México D.F.	27/11/1996	22.00
10354	Pericles Comidas clásicas, México D.F.	14/11/1996	53.80
10322	Pericles Comidas clásicas, México D.F.	04/10/1996	0.40
10319	Tortuga Restaurante, México D.F.	02/10/1996	64.50
10308	Ana Trujillo Emparedados y helados, México D.F.	18/09/1996	1.61
10304	Tortuga Restaurante, México D.F.	12/09/1996	63.79
10293	Tortuga Restaurante, México D.F.	29/08/1996	21.18
10276	Tortuga Restaurante, México D.F.	08/08/1996	13.84
10259	Centro comercial Moctezuma, México D.F.	18/07/1996	3.25
Count : 10		Total :	327.86
Count : 10		Total :	327.86

Order	Name, City	Date	Freight
<i>Countries Starting With Letter : N</i>			
Norway		Wed, Dec 18, 1996	
10387	Santé Gourmet, Stavern	18/12/1996	93.63
Count : 1		Total :	93.63
Count : 1		Total :	93.63

Count :	10	Total :	327.86
---------	----	---------	--------

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

Count :	11	5/8	Total :	421.49
---------	----	-----	---------	--------

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
-------	------------	------	---------

Countries Starting With Letter : **P**

Poland		Thu, Dec 5, 1996	
10374	Wolski Zajazd, Warszawa	05/12/1996	3.94
Count : 1		Total : 3.94	
Portugal		Mon, Mar 31, 1997	
10491	Furia Bacalhau e Frutos do Mar, Lisboa	31/03/1997	16.96
10477	Princesa Isabel Vinhos, Lisboa	17/03/1997	13.02
10464	Furia Bacalhau e Frutos do Mar, Lisboa	04/03/1997	89.00
10433	Princesa Isabel Vinhos, Lisboa	03/02/1997	73.83
10397	Princesa Isabel Vinhos, Lisboa	27/12/1996	60.26
10352	Furia Bacalhau e Frutos do Mar, Lisboa	12/11/1996	1.30
10336	Princesa Isabel Vinhos, Lisboa	23/10/1996	15.51
10328	Furia Bacalhau e Frutos do Mar, Lisboa	14/10/1996	87.03
Count : 8		Total : 356.91	
Count : 9		Total : 360.85	

Order	Name, City	Date	Freight
-------	------------	------	---------

Countries Starting With Letter : **S**

Spain		Mon, Jan 27, 1997	
10426	Galería del gastronómo, Barcelona	27/01/1997	18.69
10366	Galería del gastronómo, Barcelona	28/11/1996	10.14
10326	Bólide Comidas preparadas, Madrid	10/10/1996	77.92
10306	Romero y tomillo, Madrid	16/09/1996	7.56
10303	Godos Cocina Típica, Sevilla	11/09/1996	107.83
10282	Romero y tomillo, Madrid	15/08/1996	12.69
10281	Romero y tomillo, Madrid	14/08/1996	2.94
Count : 7		Total : 237.77	
Sweden		Fri, Feb 28, 1997	
10460	Folk och få HB, Bräcke	28/02/1997	16.27
10445	Berglunds snabbköp, Luleå	13/02/1997	9.30
10444	Berglunds snabbköp, Luleå	12/02/1997	3.50
10434	Folk och få HB, Bräcke	03/02/1997	17.92
10384	Berglunds snabbköp, Luleå	16/12/1996	168.64
10378	Folk och få HB, Bräcke	10/12/1996	5.44
10327	Folk och få HB, Bräcke	11/10/1996	63.36
10280	Berglunds snabbköp, Luleå	14/08/1996	8.98
10278	Berglunds snabbköp, Luleå	12/08/1996	92.69
10264	Folk och få HB, Bräcke	24/07/1996	3.67
Count : 10		Total : 389.77	
Switzerland		Mon, Jan 20, 1997	
10419	Richter Supermarkt, Genève	20/01/1997	137.35
10370	Chop-suey Chinese, Bern	03/12/1996	1.17
10255	Richter Supermarkt, Genève	12/07/1996	148.33
10254	Chop-suey Chinese, Bern	11/07/1996	22.98
Count : 4		Total : 309.83	
Count : 21		Total : 937.37	

Count :	9	Total :	360.85
---------	---	---------	--------

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

Count :	30	6/8	Total :	1298.22
---------	----	-----	---------	---------

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
Countries Starting With Letter : U			
UK Mon, Mar 24, 1997			
10484	B's Beverages, London	24/03/1997	6.88
10473	Island Trading, Cowes Isle of	13/03/1997	16.37
10472	Seven Seas Imports, London	12/03/1997	4.20
10471	B's Beverages, London	11/03/1997	45.59
10462	Consolidated Holdings, London	03/03/1997	6.17
10453	Around the Horn, Colchester Essex	21/02/1997	25.36
10435	Consolidated Holdings, London	04/02/1997	9.21
10400	Eastern Connection, London	01/01/1997	83.93
10388	Seven Seas Imports, London	19/12/1996	34.86
10383	Around the Horn, Colchester Essex	16/12/1996	34.24
10377	Seven Seas Imports, London	09/12/1996	22.21
10364	Eastern Connection, London	26/11/1996	71.97
10359	Seven Seas Imports, London	21/11/1996	288.43
10355	Around the Horn, Colchester Essex	15/11/1996	41.95
10321	Island Trading, Cowes Isle of	03/10/1996	3.43
10318	Island Trading, Cowes Isle of	01/10/1996	4.73
10315	Island Trading, Cowes Isle of	26/09/1996	41.76
10289	B's Beverages, London	26/08/1996	22.77
Count :	18	Total :	764.06
USA Mon, Mar 24, 1997			
10483	White Clover Markets, Seattle WA	24/03/1997	15.28
10482	Lazy K Kountry Store, Walla Walla WA	21/03/1997	7.48
10479	Rattlesnake Canyon Grocery, Albuquerque NM	19/03/1997	708.95
10469	White Clover Markets, Seattle WA	10/03/1997	60.18
10452	Save-a-lot Markets, Boise ID	20/02/1997	140.26
10441	Old World Delicatessen, Anchorage AK	10/02/1997	73.02
10440	Save-a-lot Markets, Boise ID	10/02/1997	86.53
10432	Split Rail Beer & Ale, Lander WY	31/01/1997	4.34
10415	Hungry Coyote Import Store, Elgin OR	15/01/1997	0.20
10401	Rattlesnake Canyon Grocery, Albuquerque NM	01/01/1997	12.51
10398	Save-a-lot Markets, Boise ID	30/12/1996	89.16
10394	Hungry Coyote Import Store, Elgin OR	25/12/1996	30.34
10393	Save-a-lot Markets, Boise ID	25/12/1996	126.56
10385	Split Rail Beer & Ale, Lander WY	17/12/1996	30.96
10375	Hungry Coyote Import Store, Elgin OR	06/12/1996	20.12
10369	Split Rail Beer & Ale, Lander WY	02/12/1996	195.68
10349	Split Rail Beer & Ale, Lander WY	08/11/1996	8.63
10346	Rattlesnake Canyon Grocery, NM	05/11/1996	142.08
Count :	36	Total :	2516.34

Order	Name, City	Date	Freight
Albuquerque			
10344	White Clover Markets, Seattle WA	01/11/1996	23.29
10338	Old World Delicatessen, Anchorage AK	25/10/1996	84.21
10329	Split Rail Beer & Ale, Lander WY	15/10/1996	191.67
10324	Save-a-lot Markets, Boise ID	08/10/1996	214.27
10317	Lonesome Pine Restaurant, Portland OR	30/09/1996	12.69
10316	Rattlesnake Canyon Grocery, Albuquerque NM	27/09/1996	150.15
10314	Rattlesnake Canyon Grocery, Albuquerque NM	25/09/1996	74.16
10310	The Big Cheese, Portland OR	20/09/1996	17.52
10307	Lonesome Pine Restaurant, Portland OR	17/09/1996	0.56
10305	Old World Delicatessen, Anchorage AK	13/09/1996	257.62
10294	Rattlesnake Canyon Grocery, Albuquerque NM	30/08/1996	147.26
10272	Rattlesnake Canyon Grocery, Albuquerque NM	02/08/1996	98.03
10271	Split Rail Beer & Ale, Lander WY	01/08/1996	4.54
10269	White Clover Markets, Seattle WA	31/07/1996	4.56
10262	Rattlesnake Canyon Grocery, Albuquerque NM	22/07/1996	48.29
Count :	33	Total :	3081.10
Count :	51	Total :	3845.16
Count :	15	Total :	1328.82

Count : 51

7/8

Total : 3845.16

Northwind Order List

Max order ID is : 10500

Order	Name, City	Date	Freight
-------	------------	------	---------

Countries Starting With Letter : **V**

Venezuela

Tue, Apr 8, 1997

10499	LILA-Supermercado, Barquisimeto	Lara	08/04/1997	102.02
10498	HILARION-Abastos, San Cristóbal	Táchir	07/04/1997	29.75
10490	HILARION-Abastos, San Cristóbal	Táchir	31/03/1997	210.19
10486	HILARION-Abastos, San Cristóbal	Táchir	26/03/1997	30.53
10485	LINO-Delicateses, I. de Margarita	Nueva	25/03/1997	64.45
10476	HILARION-Abastos, San Cristóbal	Táchir	17/03/1997	4.41
10461	LILA-Supermercado, Barquisimeto	Lara	28/02/1997	148.61
10405	LINO-Delicateses, I. de Margarita	Nueva	06/01/1997	34.82
10395	HILARION-Abastos, San Cristóbal	Táchir	26/12/1996	184.41
10381	LILA-Supermercado, Barquisimeto	Lara	12/12/1996	7.99
10357	LILA-Supermercado, Barquisimeto	Lara	19/11/1996	34.88
10330	LILA-Supermercado, Barquisimeto	Lara	16/10/1996	12.75
10296	LILA-Supermercado, Barquisimeto	Lara	03/09/1996	0.12
10283	LILA-Supermercado, Barquisimeto	Lara	16/08/1996	84.81
10268	GROSELLA-Restaurante, Caracas	DF	30/07/1996	66.29
10257	HILARION-Abastos, San Cristóbal	Táchir	16/07/1996	81.91
Count : 16		Total :		1097.94
Count : 16		Total :		1097.94



T h a t ' s A l l F o l k s !
Hei_remind_me_to_put myself up for
abduction. END!

Count :	16	Total :	1097.94
---------	----	---------	---------

Count : 16 8/8 Total : 1097.94