

Shipments to Argentina

		1997						1998					Total
		Jan	Feb	May	Oct	Dec	1997 Total	Jan	Feb	Mar	Apr	1998 Total	
No region	Buenos Aires	29.8	38.8	25.3	22.6	1.1	117.7	19.8	93.7	146.1	217.9	477.4	595.1
	Total	29.8	38.8	25.3	22.6	1.1	117.7	19.8	93.7	146.1	217.9	477.4	595.1
Argentina Total		29.8	38.8	25.3	22.6	1.1	117.7	19.8	93.7	146.1	217.9	477.4	595.1

Shipments to Austria

		1996				1997												1998					Total
		Jul	Nov	Dec	1996 Total	Jan	Feb	Mar	Apr	May	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar	Apr	1998 Total	
No region	Graz	286.6	162.3	323.1	772.0	141.7	506.7	0.0	0.0	790.0	122.8	477.9	78.1	272.5	0.0	593.4	2983.1	549.7	263.0	353.1	946.5	2112.2	5867.3
	Salzburg	0.0	360.6	0.0	360.6	122.5	0.0	31.3	5.3	339.2	35.1	0.0	0.0	96.5	117.3	0.0	747.2	25.2	0.0	0.0	53.0	78.3	1186.1
	Total	286.6	523.0	323.1	1132.6	264.1	506.7	31.3	5.3	1129.2	158.0	477.9	78.1	369.0	117.3	593.4	3730.3	574.9	263.0	353.1	999.5	2190.5	7053.4
Austria Total		286.6	523.0	323.1	1132.6	264.1	506.7	31.3	5.3	1129.2	158.0	477.9	78.1	369.0	117.3	593.4	3730.3	574.9	263.0	353.1	999.5	2190.5	7053.4

Shipments to Belgium

		1996			1997						1998					Total
		Jul	Oct	1996 Total	Mar	Apr	May	Aug	Dec	1997 Total	Jan	Feb	Mar	Apr	1998 Total	
No region	Bruxelles	0.0	0.0	0.0	0.0	0.0	66.7	6.2	155.6	228.5	0.0	152.7	0.0	77.7	230.4	458.9
	Charleroi	51.3	6.3	57.6	161.8	68.5	0.0	0.0	1.6	231.9	480.8	5.6	15.6	29.8	531.7	821.2
	Total	51.3	6.3	57.6	161.8	68.5	66.7	6.2	157.2	460.5	480.8	158.4	15.6	107.4	762.1	1280.1

Belgium Total	51.3	6.3	57.6	161.8	68.5	66.7	6.2	157.2	460.5	480.8	158.4	15.6	107.4	762.1	1280.1
---------------	------	-----	------	-------	------	------	-----	-------	-------	-------	-------	------	-------	-------	--------

Shipments to Brazil

		1996						1997												1998				
		Jul	Aug	Sep	Nov	Dec	1996 Total	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar	Apr
RJ	Rio de Janeiro	127.1	12.8	36.2	0.0	45.0	221.0	99.2	0.0	133.0	0.0	68.7	60.4	62.5	51.0	72.2	15.8	9.5	130.3	702.6	68.9	198.6	106.7	228.1
	Total	127.1	12.8	36.2	0.0	45.0	221.0	99.2	0.0	133.0	0.0	68.7	60.4	62.5	51.0	72.2	15.8	9.5	130.3	702.6	68.9	198.6	106.7	228.1
SP	Campinas	0.0	0.0	0.0	0.0	0.0	0.0	0.0	24.5	0.0	0.0	0.0	0.0	0.0	0.0	7.1	33.8	212.4	28.2	306.1	3.0	0.0	5.0	0.0
	Resende	14.0	0.0	0.0	0.0	0.0	14.0	44.1	0.0	0.0	0.0	0.0	0.0	13.4	0.0	0.1	0.0	0.0	0.0	57.7	60.1	0.0	63.0	0.0
	Sao Paulo	0.0	0.0	81.0	3.1	904.8	988.9	129.5	0.0	83.0	116.3	0.0	0.0	82.4	201.3	282.6	0.0	73.9	110.9	1079.9	147.0	191.3	158.9	0.0
	Total	14.0	0.0	81.0	3.1	904.8	1002.9	173.6	24.5	83.0	116.3	0.0	0.0	95.8	201.3	289.9	33.8	286.4	139.1	1443.7	210.2	191.3	226.9	0.0
Brazil Total		141.0	12.8	117.2	3.1	949.8	1223.9	272.9	24.5	216.0	116.3	68.7	60.4	158.3	252.3	362.1	49.5	295.9	269.4	2146.3	279.0	389.9	333.6	228.1

		1998		Total
		May	1998 Total	
RJ	Rio de Janeiro	73.5	675.8	1599.5
	Total	73.5	675.8	1599.5
SP	Campinas	8.3	16.3	322.4
	Resende	0.0	123.1	194.7

SP	Sao Paulo	30.0	527.2	2596.1
	Total	38.3	666.6	3113.2
Brazil Total		111.9	1342.5	4712.6

Shipments to Canada

		1996				1997								1998				Total
		Oct	Nov	Dec	1996 Total	Jan	Feb	Apr	Jun	Jul	Aug	Nov	1997 Total	Jan	Mar	Apr	1998 Total	
BC	Tsawassen	0.0	0.0	47.4	47.4	26.0	44.2	62.9	0.0	0.0	0.0	243.7	376.8	0.0	208.5	90.7	299.1	723.4
	Vancouver	0.0	0.0	0.0	0.0	0.0	0.0	4.7	0.0	0.0	0.9	0.0	5.6	4.3	0.0	0.0	4.3	9.9
	Total	0.0	0.0	47.4	47.4	26.0	44.2	67.5	0.0	0.0	0.9	243.7	382.4	4.3	208.5	90.7	303.4	733.3
Québec	Montréal	52.8	15.7	20.4	88.9	370.6	4.1	7.1	196.1	423.9	245.7	57.8	1305.3	0.0	0.0	0.0	0.0	1394.2
	Total	52.8	15.7	20.4	88.9	370.6	4.1	7.1	196.1	423.9	245.7	57.8	1305.3	0.0	0.0	0.0	0.0	1394.2
Canada Total		52.8	15.7	67.8	136.3	396.7	48.2	74.7	196.1	423.9	246.7	301.5	1687.8	4.3	208.5	90.7	303.4	2127.5

Shipments to Denmark

		1996			1997								1998				Total	
		Nov	Dec	1996 Total	Jan	Mar	Jun	Jul	Sep	Oct	Nov	Dec	1997 Total	Jan	Mar	Apr		1998 Total
No region	Kobenhavn	26.8	0.0	26.8	70.3	0.0	9.8	0.0	66.3	0.0	0.0	0.0	146.4	257.3	0.0	0.0	257.3	430.4
	Århus	0.0	13.6	13.6	27.4	145.0	0.0	58.8	0.0	299.1	69.2	65.1	664.6	0.0	203.7	65.5	269.2	947.3
	Total	26.8	13.6	40.3	97.7	145.0	9.8	58.8	66.3	299.1	69.2	65.1	810.9	257.3	203.7	65.5	526.5	1377.8
Denmark Total		26.8	13.6	40.3	97.7	145.0	9.8	58.8	66.3	299.1	69.2	65.1	810.9	257.3	203.7	65.5	526.5	1377.8

Shipments to Finland

		1996				1997												1998				Total
		Jul	Aug	Oct	1996 Total	Jan	Feb	Mar	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Feb	Mar	Apr	1998 Total	
No region	Helsinki	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.8	22.8	16.7	0.0	0.0	40.2	9.3	38.1	0.8	48.2	88.4
	Oulu	25.7	136.5	35.2	197.4	26.5	20.0	180.4	58.6	149.5	7.3	1.1	0.0	0.0	79.3	73.2	595.9	0.0	0.0	29.2	29.2	822.5
	Total	25.7	136.5	35.2	197.4	26.5	20.0	180.4	58.6	149.5	7.3	1.9	22.8	16.7	79.3	73.2	636.1	9.3	38.1	29.9	77.3	910.9
Finland Total		25.7	136.5	35.2	197.4	26.5	20.0	180.4	58.6	149.5	7.3	1.9	22.8	16.7	79.3	73.2	636.1	9.3	38.1	29.9	77.3	910.9

Shipments to France

		1996							1997													1998		
		Jul	Aug	Sep	Oct	Nov	Dec	1996 Total	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar
No region	Lille	0.0	0.0	0.0	0.0	0.0	0.0	0.0	11.3	0.0	1.4	0.0	0.0	0.0	0.0	487.4	0.0	0.0	0.0	137.9	637.9	0.0	0.0	0.0
	Lyon	41.3	0.0	0.0	8.6	0.0	0.0	49.9	0.0	25.1	12.0	0.0	194.7	0.0	0.0	0.0	0.0	0.0	0.0	0.0	231.9	211.5	0.0	0.0
	Marseille	0.0	0.0	0.0	10.2	262.4	0.0	272.5	0.0	0.0	64.6	350.6	11.1	0.0	0.0	0.0	0.0	176.4	53.8	0.0	656.4	0.0	236.2	154.4
	Nantes	0.0	0.0	24.7	0.0	0.0	0.0	24.7	0.0	0.0	0.0	0.0	0.0	0.0	1.9	0.0	30.3	4.4	0.0	0.0	36.6	0.0	52.0	0.0
	Paris	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	2.9	0.0	2.9	0.0	9.2	87.4
	Reims	32.4	6.0	1.1	0.0	0.0	0.0	39.5	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	18.9	0.0	18.9	0.0	0.0	0.0
	Strasbourg	0.0	55.3	5.7	0.0	0.0	131.7	192.7	0.0	210.0	0.0	0.0	0.0	96.5	59.1	30.4	27.9	0.0	0.0	0.0	423.8	0.0	7.1	0.0
	Toulouse	0.0	0.0	0.0	0.0	19.6	64.6	84.3	95.7	10.7	0.0	53.3	0.0	0.0	0.0	27.7	0.0	0.0	0.0	249.9	437.2	43.3	0.0	68.3
	Versailles	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	52.5	15.2
	Total	73.7	61.3	31.6	18.8	282.0	196.3	663.7	106.9	245.7	77.9	404.0	205.8	96.5	61.0	545.4	58.3	180.8	75.6	387.9	2445.6	254.8	357.0	325.2

France Total		73.7	61.3	31.6	18.8	282.0	196.3	663.7	106.9	245.7	77.9	404.0	205.8	96.5	61.0	545.4	58.3	180.8	75.6	387.9	2445.6	254.8	357.0	325.2
		1998		Total																				
		Apr	1998 Total																					
No region	Lille	0.0	0.0	637.9																				
	Lyon	0.0	211.5	493.2																				
	Marseille	0.0	390.6	1319.6																				
	Nantes	121.8	173.8	235.1																				
	Paris	8.8	105.4	108.3																				
	Reims	0.0	0.0	58.4																				
	Strasbourg	0.0	7.1	623.7																				
	Toulouse	0.0	111.5	633.0																				
	Versailles	19.8	87.5	87.5																				
	Total	150.4	1087.5	4196.8																				
France Total		150.4	1087.5	4196.8																				

Shipments to Germany

		1996							1997												1998			
		Jul	Aug	Sep	Oct	Nov	Dec	1996 Total	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar
No region	Aachen	0.0	0.0	0.0	0.0	0.0	36.0	36.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	112.6	0.0	0.0
	Berlin	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	29.5	85.0	0.0	0.0	114.4	69.5	0.0	40.4
	Brandenburg	0.0	0.0	0.0	69.7	0.0	0.0	69.7	0.0	8.1	55.7	0.0	32.1	0.0	0.0	32.3	0.0	170.9	0.0	0.0	299.2	337.4	77.8	0.0

No region	Cunewalde	0.0	382.1	0.0	2.0	249.1	183.2	816.3	17.5	0.0	189.1	0.0	417.6	1007.6	194.7	0.0	364.1	1257.3	3.5	42.7	3494.3	255.7	394.8	307.7
	Frankfurt a.M.	0.0	102.4	0.0	0.0	110.4	0.0	212.8	0.0	0.0	0.0	36.2	73.3	155.6	32.1	174.2	0.0	0.0	0.0	91.3	562.7	0.0	73.6	32.0
	Köln	55.1	0.0	0.0	0.0	0.0	0.0	55.1	91.5	0.0	0.0	0.0	5.0	121.0	75.9	0.0	145.6	0.0	0.0	157.6	596.5	71.5	0.0	0.0
	Leipzig	0.0	125.8	0.0	0.0	0.0	0.0	125.8	0.0	0.0	0.0	0.0	0.0	127.3	0.0	0.0	0.0	0.6	0.0	0.0	127.9	58.1	0.0	10.2
	Mannheim	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	9.0	0.0	0.0	27.7	1.9	0.0	0.0	0.0	0.0	38.6	0.0	53.8	44.7
	München	0.0	208.6	0.0	108.3	54.8	0.0	371.7	135.4	0.0	0.0	4.9	0.0	36.7	0.0	97.2	328.6	59.2	0.0	0.0	661.9	16.9	76.1	33.9
	Münster	11.6	0.0	0.0	0.0	0.0	0.0	11.6	0.0	22.9	0.0	0.0	0.0	1.4	0.0	27.8	0.0	0.0	0.0	0.0	52.1	0.0	0.0	0.0
	Stuttgart	0.0	0.0	45.1	40.3	37.5	0.0	122.8	0.0	0.0	0.0	105.7	0.0	0.0	0.0	64.9	67.8	0.0	0.0	0.0	238.4	0.0	0.0	0.0
	Total	66.7	818.9	45.1	220.2	451.8	219.2	1821.8	244.4	31.0	244.8	155.8	528.0	1449.6	330.4	398.4	935.6	1573.0	3.5	291.6	6186.1	921.7	676.1	468.9
Germany Total		66.7	818.9	45.1	220.2	451.8	219.2	1821.8	244.4	31.0	244.8	155.8	528.0	1449.6	330.4	398.4	935.6	1573.0	3.5	291.6	6186.1	921.7	676.1	468.9

		1998			Total
		Apr	May	1998 Total	
No region	Aachen	149.5	8.0	270.1	306.0
	Berlin	1.2	0.0	111.2	225.6
	Brandenburg	29.6	0.0	444.8	813.7
	Cunewalde	336.8	0.0	1295.0	5605.6
	Frankfurt a.M.	0.0	0.0	105.6	881.0
	Köln	139.6	0.0	211.1	862.7
	Leipzig	0.0	0.0	68.3	322.0
	Mannheim	0.0	0.0	98.5	137.1

No region	München	242.9	0.0	369.8	1403.4
	Münster	62.2	0.0	62.2	126.0
	Stuttgart	71.6	0.0	71.6	432.9
	Total	1033.5	8.0	3108.2	11116.1
Germany Total		1033.5	8.0	3108.2	11116.1

Shipments to Ireland

		1996				1997									1998					Total
		Sep	Oct	Dec	1996 Total	Jan	Feb	Apr	May	Jun	Sep	Oct	Nov	1997 Total	Feb	Mar	Apr	May	1998 Total	
Co. Cork	Cork	168.2	89.4	124.1	381.8	35.0	56.6	16.7	62.8	34.0	159.9	606.7	44.1	1015.8	603.5	580.9	91.5	81.7	1357.7	2755.2
	Total	168.2	89.4	124.1	381.8	35.0	56.6	16.7	62.8	34.0	159.9	606.7	44.1	1015.8	603.5	580.9	91.5	81.7	1357.7	2755.2
Ireland Total		168.2	89.4	124.1	381.8	35.0	56.6	16.7	62.8	34.0	159.9	606.7	44.1	1015.8	603.5	580.9	91.5	81.7	1357.7	2755.2

Shipments to Italy

		1996			1997										1998					Total	
		Aug	Sep	1996 Total	Jan	Feb	Mar	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Mar	Apr	May		1998 Total
No region	Bergamo	26.9	17.7	44.6	156.0	0.0	4.9	0.0	0.0	47.5	0.0	0.0	2.4	70.1	280.8	65.5	78.8	0.0	0.0	144.3	469.8
	Reggio Emilia	0.0	7.4	7.4	0.0	25.0	0.0	23.0	0.5	0.0	4.4	0.0	0.0	89.9	142.8	59.8	50.9	28.7	0.0	139.4	289.6
	Torino	0.0	0.0	0.0	3.0	0.0	0.0	0.0	0.0	0.0	0.0	5.0	7.7	0.0	15.7	1.4	0.0	47.1	11.0	59.4	75.1
	Total	26.9	25.1	52.1	159.0	25.0	4.9	23.0	0.5	47.5	4.4	5.0	10.1	160.0	439.3	126.6	129.7	75.8	11.0	343.1	834.5
Italy Total		26.9	25.1	52.1	159.0	25.0	4.9	23.0	0.5	47.5	4.4	5.0	10.1	160.0	439.3	126.6	129.7	75.8	11.0	343.1	834.5

Shipments to Mexico

		1996							1997								1998						
		Jul	Aug	Sep	Oct	Nov	Dec	1996 Total	Mar	Apr	May	Jun	Aug	Sep	Oct	Dec	1997 Total	Jan	Feb	Mar	Apr	May	1998 Total
No region	México D.F.	3.2	13.8	86.6	64.9	53.8	22.0	244.4	83.5	116.8	233.8	103.4	43.9	6.0	36.1	12.0	635.5	54.4	58.4	43.4	46.0	15.7	217.9
	Total	3.2	13.8	86.6	64.9	53.8	22.0	244.4	83.5	116.8	233.8	103.4	43.9	6.0	36.1	12.0	635.5	54.4	58.4	43.4	46.0	15.7	217.9
Mexico Total		3.2	13.8	86.6	64.9	53.8	22.0	244.4	83.5	116.8	233.8	103.4	43.9	6.0	36.1	12.0	635.5	54.4	58.4	43.4	46.0	15.7	217.9

Total
-------

No region	México D.F.	1097.8
	Total	1097.8
Mexico Total		1097.8

Shipments to Norway

		1996		1997			1998				Total
		Dec	1996 Total	May	Aug	1997 Total	Jan	Mar	Apr	1998 Total	
No region	Stavern	93.6	93.6	13.4	38.6	52.0	72.2	53.0	4.6	129.9	275.5
	Total	93.6	93.6	13.4	38.6	52.0	72.2	53.0	4.6	129.9	275.5
Norway Total		93.6	93.6	13.4	38.6	52.0	72.2	53.0	4.6	129.9	275.5



Shipments to Poland

		1996		1997			1998					Total
		Dec	1996 Total	Aug	Dec	1997 Total	Feb	Mar	Apr	May	1998 Total	
No region	Warszawa	3.9	3.9	80.7	23.8	104.4	12.0	26.3	20.3	8.7	67.4	175.7
	Total	3.9	3.9	80.7	23.8	104.4	12.0	26.3	20.3	8.7	67.4	175.7
Poland Total		3.9	3.9	80.7	23.8	104.4	12.0	26.3	20.3	8.7	67.4	175.7

Shipments to Portugal

		1996			1997							1998			Total
		Oct	Nov	1996 Total	Jan	Mar	Apr	Jun	Jul	Sep	1997 Total	Mar	Apr	1998 Total	
No region	Lisboa	102.5	1.3	103.8	60.3	175.9	17.0	72.9	7.5	1.3	334.8	2.7	202.2	204.9	643.5
	Total	102.5	1.3	103.8	60.3	175.9	17.0	72.9	7.5	1.3	334.8	2.7	202.2	204.9	643.5
Portugal Total		102.5	1.3	103.8	60.3	175.9	17.0	72.9	7.5	1.3	334.8	2.7	202.2	204.9	643.5

Shipments to Spain

		1996					1997						1998				Total
		Aug	Sep	Oct	Dec	1996 Total	Feb	Jun	Jul	Aug	Dec	1997 Total	Feb	Mar	Apr	1998 Total	
No region	Barcelona	0.0	0.0	0.0	10.1	10.1	18.7	0.0	6.5	0.0	0.0	25.2	1.2	1.4	0.0	2.6	38.0
	Madrid	15.6	7.6	77.9	0.0	101.1	0.0	0.0	0.0	0.0	97.1	97.1	0.0	8.3	49.2	57.4	255.6
	Sevilla	0.0	107.8	0.0	0.0	107.8	0.0	4.3	0.0	85.5	0.0	89.8	246.8	61.6	62.3	370.7	568.3
	Total	15.6	115.4	77.9	10.1	219.1	18.7	4.3	6.5	85.5	97.1	212.1	248.0	71.2	111.5	430.7	861.9

Spain Total	15.6	115.4	77.9	10.1	219.1	18.7	4.3	6.5	85.5	97.1	212.1	248.0	71.2	111.5	430.7	861.9
-------------	------	-------	------	------	-------	------	-----	-----	------	------	-------	-------	------	-------	-------	-------

Shipments to Sweden

		1996					1997										1998						Total
		Aug	Sep	Oct	Dec	1996 Total	Feb	Mar	May	Jun	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar	Apr	May	1998 Total	
No region	Bräcke	3.7	0.0	63.4	5.4	72.5	17.9	16.3	188.0	242.2	0.0	0.0	152.3	0.0	376.9	993.7	1.2	88.0	47.4	415.9	59.4	611.9	1678.1
	Luleå	92.7	9.0	0.0	168.6	270.3	12.8	0.0	244.8	116.4	138.7	151.0	13.4	110.1	6.8	794.0	13.3	298.0	32.4	151.5	0.0	495.2	1559.5
	Total	96.4	9.0	63.4	174.1	342.8	30.7	16.3	432.8	358.6	138.7	151.0	165.7	110.1	383.7	1787.7	14.5	386.0	79.8	567.4	59.4	1107.1	3237.6
Sweden Total		96.4	9.0	63.4	174.1	342.8	30.7	16.3	432.8	358.6	138.7	151.0	165.7	110.1	383.7	1787.7	14.5	386.0	79.8	567.4	59.4	1107.1	3237.6

Shipments to Switzerland

		1996			1997						1998			Total
		Jul	Dec	1996 Total	Jan	May	Sep	Nov	Dec	1997 Total	Mar	Apr	1998 Total	
No region	Bern	23.0	1.2	24.1	0.0	91.8	0.0	128.1	0.0	219.8	0.0	123.2	123.2	367.2
	Genève	148.3	0.0	148.3	137.4	78.8	232.4	0.0	269.0	717.6	13.6	115.6	129.2	995.1
	Total	171.3	1.2	172.5	137.4	170.6	232.4	128.1	269.0	937.4	13.6	238.8	252.4	1362.3
Switzerland Total		171.3	1.2	172.5	137.4	170.6	232.4	128.1	269.0	937.4	13.6	238.8	252.4	1362.3

Shipments to UK

		1996					1997													1998			
		Aug	Oct	Nov	Dec	1996 Total	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar	Apr
Essex	Colchester	0.0	0.0	42.0	34.2	76.2	0.0	25.4	0.0	0.0	0.0	73.0	0.0	0.0	0.0	21.7	34.7	146.3	301.1	4.5	3.0	53.3	33.8
	Total	0.0	0.0	42.0	34.2	76.2	0.0	25.4	0.0	0.0	0.0	73.0	0.0	0.0	0.0	21.7	34.7	146.3	301.1	4.5	3.0	53.3	33.8
Isle of Wight	Cowes	0.0	49.9	0.0	0.0	49.9	0.0	0.0	16.4	0.0	0.0	0.0	0.0	23.7	0.9	0.0	0.0	61.5	102.5	157.1	0.0	54.2	0.0
	Total	0.0	49.9	0.0	0.0	49.9	0.0	0.0	16.4	0.0	0.0	0.0	0.0	23.7	0.9	0.0	0.0	61.5	102.5	157.1	0.0	54.2	0.0
No region	London	22.8	0.0	288.4	129.0	440.2	83.9	9.2	56.0	39.0	169.3	178.4	59.6	0.0	0.0	0.0	1.4	16.6	613.3	203.0	143.3	5.4	383.7
	Total	22.8	0.0	288.4	129.0	440.2	83.9	9.2	56.0	39.0	169.3	178.4	59.6	0.0	0.0	0.0	1.4	16.6	613.3	203.0	143.3	5.4	383.7
UK Total		22.8	49.9	330.4	163.3	566.3	83.9	34.6	72.3	39.0	169.3	251.4	59.6	23.7	0.9	21.7	36.1	224.4	1016.9	364.6	146.3	112.9	417.5

		1998		Total
		May	1998 Total	
Essex	Colchester	0.0	94.7	472.0
	Total	0.0	94.7	472.0
Isle of Wight	Cowes	0.0	211.2	363.7
	Total	0.0	211.2	363.7
No region	London	329.7	1065.1	2118.7
	Total	329.7	1065.1	2118.7
UK Total		329.7	1371.0	2954.3

Shipments to USA

		1996							1997													1998		
		Jul	Aug	Sep	Oct	Nov	Dec	1996 Total	Jan	Feb	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar
AK	Anchorage	0.0	0.0	0.0	341.8	0.0	0.0	341.8	0.0	0.0	73.0	0.0	0.0	0.0	5.2	0.0	26.6	135.6	0.0	0.0	240.5	45.5	171.0	144.4
	Total	0.0	0.0	0.0	341.8	0.0	0.0	341.8	0.0	0.0	73.0	0.0	0.0	0.0	5.2	0.0	26.6	135.6	0.0	0.0	240.5	45.5	171.0	144.4
CA	San Francisco	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	13.7	0.0	0.0	0.0	97.4	0.0	111.1	0.0	91.0	0.0
	Total	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	13.7	0.0	0.0	0.0	97.4	0.0	111.1	0.0	91.0	0.0
ID	Boise	0.0	0.0	0.0	214.3	0.0	0.0	214.3	215.7	226.8	0.0	367.6	0.0	252.5	200.2	700.3	352.7	698.0	307.1	8.2	3329.2	14.6	626.8	400.8
	Total	0.0	0.0	0.0	214.3	0.0	0.0	214.3	215.7	226.8	0.0	367.6	0.0	252.5	200.2	700.3	352.7	698.0	307.1	8.2	3329.2	14.6	626.8	400.8
MT	Butte	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	94.8	0.0	0.0	0.0	20.2	115.1	0.0	0.0	0.0
	Total	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	94.8	0.0	0.0	0.0	20.2	115.1	0.0	0.0	0.0
NM	Albuquerque	48.3	98.0	147.3	224.3	142.1	0.0	660.0	12.5	0.0	709.0	0.0	0.0	13.8	103.4	0.0	0.0	0.0	0.0	18.7	857.3	211.6	280.6	0.0
	Total	48.3	98.0	147.3	224.3	142.1	0.0	660.0	12.5	0.0	709.0	0.0	0.0	13.8	103.4	0.0	0.0	0.0	0.0	18.7	857.3	211.6	280.6	0.0
OR	Elgin	0.0	0.0	0.0	0.0	0.0	20.1	20.1	30.5	0.0	0.0	0.0	0.0	0.0	45.1	0.0	0.0	111.3	0.0	0.0	187.0	0.0	0.0	0.0
	Eugene	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	3.3	0.0	4.4	135.1	133.3	0.0	0.0	0.0	276.1	0.0	719.8	33.7
	Portland	0.0	0.0	18.1	12.7	0.0	0.0	30.8	0.0	0.0	0.0	0.0	24.9	0.0	0.0	0.0	27.6	0.0	3.0	0.0	55.5	237.3	2.5	0.0
	Total	0.0	0.0	18.1	12.7	0.0	20.1	50.9	30.5	0.0	0.0	0.0	28.3	0.0	49.6	135.1	160.9	111.3	3.0	0.0	518.5	237.3	722.2	33.7
WA	Kirkland	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	63.0	0.0	0.0	0.0	0.0	0.0	0.0	63.0	7.0	0.0	0.0
	Seattle	0.0	4.6	0.0	0.0	23.3	0.0	27.9	0.0	0.0	60.2	74.4	0.0	0.0	0.0	16.3	0.0	241.9	103.6	0.0	496.4	0.0	177.9	0.0
	Walla Walla	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	7.5	0.0	11.9	0.0	0.0	0.0	0.0	0.0	0.0	19.4	0.0	0.0	0.0
	Total	0.0	4.6	0.0	0.0	23.3	0.0	27.9	0.0	0.0	60.2	81.9	0.0	74.9	0.0	16.3	0.0	241.9	103.6	0.0	578.8	7.0	177.9	0.0

WY	Lander	0.0	4.5	0.0	191.7	8.6	226.6	431.5	0.0	4.3	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	73.2	77.6	36.7	0.0	0.0
	Total	0.0	4.5	0.0	191.7	8.6	226.6	431.5	0.0	4.3	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	73.2	77.6	36.7	0.0	0.0
USA Total		48.3	107.1	165.3	984.8	174.0	246.8	1726.3	258.8	231.1	842.2	449.5	28.3	341.2	372.2	946.5	540.2	1186.8	511.1	120.3	5828.1	552.7	2069.5	578.9

1998			Total
Apr	May	1998 Total	

AK	Anchorage	40.3	0.0	401.2	983.5
	Total	40.3	0.0	401.2	983.5
CA	San Francisco	0.0	0.0	91.0	202.1
	Total	0.0	0.0	91.0	202.1
ID	Boise	2067.9	30.1	3140.2	6683.7
	Total	2067.9	30.1	3140.2	6683.7
MT	Butte	14.9	0.0	14.9	130.0
	Total	14.9	0.0	14.9	130.0
NM	Albuquerque	116.3	0.0	608.4	2125.7
	Total	116.3	0.0	608.4	2125.7
OR	Elgin	0.0	0.0	0.0	207.1
	Eugene	25.2	0.0	778.7	1054.8
	Portland	15.9	0.0	255.7	342.0
	Total	41.1	0.0	1034.4	1603.8

WA	Kirkland	0.0	0.0	7.0	70.0
	Seattle	606.2	44.7	828.8	1353.1
	Walla Walla	0.0	0.0	0.0	19.4
	Total	606.2	44.7	835.8	1442.5
WY	Lander	13.0	0.0	49.6	558.7
	Total	13.0	0.0	49.6	558.7
USA Total		2899.6	74.8	6175.5	13729.9

Shipments to Venezuela

		1996						1997												1998			
		Jul	Aug	Sep	Oct	Dec	1996 Total	Jan	Mar	Apr	May	Jun	Jul	Aug	Sep	Oct	Nov	Dec	1997 Total	Jan	Feb	Mar	Apr
DF	Caracas	0.0	66.3	0.0	0.0	0.0	66.3	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	1.5	1.5	0.0	0.0	0.0	0.0
	Total	0.0	66.3	0.0	0.0	0.0	66.3	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	0.0	1.5	1.5	0.0	0.0	0.0	0.0
Lara	Barquisimeto	0.0	84.8	0.1	12.8	42.9	140.6	0.0	148.6	102.0	48.2	0.0	0.0	0.0	0.0	0.0	0.0	42.1	340.9	164.0	1.2	0.0	73.9
	Total	0.0	84.8	0.1	12.8	42.9	140.6	0.0	148.6	102.0	48.2	0.0	0.0	0.0	0.0	0.0	0.0	42.1	340.9	164.0	1.2	0.0	73.9
Nueva Esparta	I. de Margarita	0.0	0.0	0.0	0.0	0.0	0.0	34.8	64.4	0.0	0.0	0.0	0.0	0.0	158.4	45.5	141.1	0.0	444.3	90.5	2.7	47.7	23.6
	Total	0.0	0.0	0.0	0.0	0.0	0.0	34.8	64.4	0.0	0.0	0.0	0.0	0.0	158.4	45.5	141.1	0.0	444.3	90.5	2.7	47.7	23.6
Táchira	San Cristóbal	81.9	0.0	0.0	0.0	0.0	81.9	184.4	4.4	270.5	0.0	83.2	58.3	187.7	0.0	0.0	3.5	0.0	792.1	26.5	92.3	105.4	40.1
	Total	81.9	0.0	0.0	0.0	0.0	81.9	184.4	4.4	270.5	0.0	83.2	58.3	187.7	0.0	0.0	3.5	0.0	792.1	26.5	92.3	105.4	40.1
Venezuela Total		81.9	151.1	0.1	12.8	42.9	288.8	219.2	217.5	372.5	48.2	83.2	58.3	187.7	158.4	45.5	144.6	43.6	1578.8	281.0	96.3	153.1	137.6

		1998		Total
		May	1998 Total	
DF	Caracas	0.0	0.0	67.8
	Total	0.0	0.0	67.8
Lara	Barquisimeto	0.0	239.1	720.6
	Total	0.0	239.1	720.6
Nueva Esparta	I. de Margarita	0.0	164.5	608.8
	Total	0.0	164.5	608.8
Táchira	San Cristóbal	120.9	385.2	1259.2
	Total	120.9	385.2	1259.2
Venezuela Total		120.9	788.8	2656.3