

# Master Report

## Total order value: 14243.4

### City List

#### Berne 2 addresses

Title Products ordered by people in Berne				Title Addresses in Berne	
Page Header				Page Header	
ID	Name	Qty	Price	ID	Name Street
0	Iron Iron	3	8.10	9	James Schneider 277 Seventh Av.
1	Chair Shoe	4	37.20	22	Bill Ott 250 - 20th Ave.
4	Ice Tea Shoe	48	57.60	Summary Count : 2	
5	Clock Clock	12	35.40	Column Footer	
6	Ice Tea Chair	23	14.70	Page Footer Page 1 of 1	
7	Telephone Shoe	5	12.60		
8	Ice Tea Clock	14	33.90		
10	Telephone Ice Tea	2	30.60		
11	Telephone Iron	17	13.20		
12	Clock Ice Tea	17	50.40		
14	Telephone Iron	15	18.60		
23	Shoe Chair	10	11.40		
24	Chair Shoe	16	10.80		
26	Shoe Shoe	39	75.60		
30	Shoe Iron	19	34.80		
31	Ice Tea Telephone	2	7.20		
33	Iron Chair	10	27.30		
35	Telephone Shoe	17	11.40		
36	Ice Tea Iron	24	4.80		
38	Clock Ice Tea	3	32.40		
41	Clock Ice Tea	10	30.90		
48	Clock Clock	4	31.50		
49	Iron Iron	23	3.30		
Summary				Count : 23	
Column Footer				Total : 337 593.70	
Page Footer				Page 1 of 1	

#### Boston 2 addresses

Title Products ordered by people in Boston						Title Addresses in Boston	
Page Header						Page Header	
ID	Name	Qty	Price	ID	Name	Qty	Price
0	Iron Iron	45	24.30	7	Telephone Shoe	24	12.60
1	Chair Shoe	24	37.20	9	Clock Telephone	4	25.80
3	Chair Chair	14	38.10	11	Telephone Iron	9	13.20
4	Ice Tea Shoe	18	19.20	12	Clock Ice Tea	4	50.40
6	Ice Tea Chair	48	44.10	13	Telephone Clock	12	27.00
Column Footer						Summary Count : 2	
Column Footer						Column Footer	
Page Footer						Page Footer Page 1 of 1	
Column Footer						Total : 149 162.90	
Page Footer						Total : 202 291.90	
Page Footer						Page 1 of 2	

## City List

### Boston (continued)

Page Header

ID	Name	Qty	Price
14	Telephone Iron	15	18.60
16	Ice Tea Shoe	1	29.10
17	Clock Ice Tea	33	66.00
18	Chair Clock	13	51.60
19	Ice Tea Ice Tea	37	33.00
20	Ice Tea Telephone	11	30.00
21	Chair Chair	32	34.20
22	Iron Iron	32	19.80
23	Shoe Chair	5	11.40
24	Chair Shoe	18	10.80
25	Shoe Shoe	36	72.90
28	Clock Iron	39	9.00
29	Chair Telephone	21	33.60
30	Shoe Iron	25	104.40
31	Ice Tea Telephone	30	14.40
33	Iron Chair	20	27.30
34	Chair Iron	17	38.40
36	Ice Tea Iron	45	9.60
37	Clock Shoe	17	14.10
38	Clock Ice Tea	18	32.40
39	Ice Tea Shoe	2	23.10
40	Shoe Clock	15	4.20
41	Clock Ice Tea	41	61.80
42	Iron Chair	21	29.70
43	Telephone Clock	8	14.10
46	Ice Tea Clock	19	9.30
47	Ice Tea Iron	14	26.70
48	Clock Clock	52	126.00
49	Iron Iron	28	6.60

Summary Count : 39

Column Footer Total : 867 1254.00

Page Footer

Page 2 of 2

### Chicago 3 addresses

Title Products ordered by people in Chicago

Title Addresses in Chicago

Page Header

Page Header

ID	Name	Qty	Price	ID	Name	Qty	Price	ID	Name	Street
0	Iron Iron	36	16.20	9	Clock Telephone	22	25.80	11	Julia White	412 Upland Pl.
1	Chair Shoe	45	148.80	10	Telephone Ice Tea	20	30.60	35	George Karsen	412 College Av.
4	Ice Tea Shoe	20	19.20	11	Telephone Iron	31	26.40	39	Mary Karsen	202 College Av.
5	Clock Clock	19	35.40	12	Clock Ice Tea	3	25.20	Summary		Count : 3
6	Ice Tea Chair	12	29.40	13	Telephone Clock	77	108.00	Column Footer		

Column Footer Total : 132 249.00 Column Footer Total : 285 465.00

Page Footer Page 1 of 1

Page Footer

Page 1 of 2

## City List

### Chicago (continued)

Page Header

ID	Name	Qty	Price
14	Telephone Iron	4	37.20
15	Ice Tea Chair	51	42.30
16	Ice Tea Shoe	10	29.10
17	Clock Ice Tea	15	66.00
19	Ice Tea Ice Tea	28	33.00
22	Iron Iron	42	19.80
23	Shoe Chair	7	11.40
24	Chair Shoe	58	43.20
25	Shoe Shoe	15	24.30
26	Shoe Shoe	68	151.20
27	Telephone Iron	8	34.50
28	Clock Iron	14	9.00
29	Chair Telephone	24	16.80
30	Shoe Iron	5	34.80
32	Clock Iron	35	114.00
33	Iron Chair	16	27.30
34	Chair Iron	23	76.80
35	Telephone Shoe	13	11.40
36	Ice Tea Iron	8	4.80
38	Clock Ice Tea	18	32.40
40	Shoe Clock	1	4.20
41	Clock Ice Tea	41	61.80
42	Iron Chair	36	59.40
43	Telephone Clock	4	14.10
44	Ice Tea Ice Tea	38	28.80
45	Iron Ice Tea	41	54.00
46	Ice Tea Clock	9	9.30
47	Ice Tea Iron	16	26.70
48	Clock Clock	11	31.50
49	Iron Iron	4	3.30

Summary Count : 40

Column Footer Total : 948 1577.40

Page Footer Page 2 of2

### Dallas 9 addresses

Title Products ordered by people in Dallas

Page Header

ID	Name	Qty	Price	ID	Name	Qty	Price
0	Iron Iron	24	16.20	4	Ice Tea Shoe	119	153.60
1	Chair Shoe	44	111.60	5	Clock Clock	56	141.60
2	Telephone Clock	2	37.20	6	Ice Tea Chair	36	44.10
3	Chair Chair	26	114.30	7	Telephone Shoe	68	50.40

Column Footer Total : 96 279.30 Column Footer Total : 375 669.00

Page Footer Page 1 of2

Title Addresses in Dallas

Page Header

ID	Name	Street
0	Laura Steel	429 Seventh Av.
4	Sylvia Ringer	365 College Av.
10	Anne Fuller	135 Upland Pl.
19	Susanne Heiniger	86 - 20th Ave.

Column Footer

Page Footer Page 1 of2

## City List

### Dallas (continued)

Page Header

ID	Name	Qty	Price
8	Ice Tea Clock	37	135.60
9	Clock Telephone	58	103.20
10	Telephone Ice Tea	22	61.20
11	Telephone Iron	4	13.20
12	Clock Ice Tea	47	100.80
13	Telephone Clock	17	27.00
14	Telephone Iron	89	111.60
15	Ice Tea Chair	48	56.40
16	Ice Tea Shoe	68	145.50
17	Clock Ice Tea	38	99.00
18	Chair Clock	13	25.80
19	Ice Tea Ice Tea	28	66.00
20	Ice Tea Telephone	37	90.00
21	Chair Chair	6	17.10
22	Iron Iron	50	39.60
23	Shoe Chair	28	34.20
24	Chair Shoe	57	54.00
25	Shoe Shoe	41	72.90
26	Shoe Shoe	49	151.20
27	Telephone Iron	28	103.50
28	Clock Iron	38	13.50
29	Chair Telephone	35	67.20
30	Shoe Iron	34	139.20
31	Ice Tea Telephone	33	21.60
32	Clock Iron	33	85.50
33	Iron Chair	46	109.20
34	Chair Iron	18	38.40
35	Telephone Shoe	17	22.80
36	Ice Tea Iron	23	14.40
37	Clock Shoe	34	28.20
38	Clock Ice Tea	28	64.80
39	Ice Tea Shoe	3	23.10
40	Shoe Clock	48	12.60
41	Clock Ice Tea	27	61.80
43	Telephone Clock	10	14.10
44	Ice Tea Ice Tea	63	86.40
45	Iron Ice Tea	24	27.00
46	Ice Tea Clock	24	9.30
47	Ice Tea Iron	38	106.80
48	Clock Clock	38	94.50

Summary Count : 48

Column Footer Total : 1754 3217.20

Page Footer

Page 2 of 2

Page Header

ID	Name	Street
36	John Steel	276 Upland Pl.
37	Michael Clancy	19 Seventh Av.
40	Susanne Miller	440 - 20th Ave.
43	Susanne Smith	2 Upland Pl.
47	Janet Fuller	445 Upland Pl.

Summary Count : 9

Column Footer

Page Footer

Page 2 of 2

## City List

### Lyon

4 addresses

Products ordered by people in Lyon				Addresses in Lyon	
Title				Title	
Page Header				Page Header	
ID	Name	Qty	Price	ID	Name
2	Telephone Clock	22	37.20	2	Anne Miller
3	Chair Chair	12	38.10	17	Laura Ott
4	Ice Tea Shoe	14	19.20	28	Susanne White
6	Ice Tea Chair	2	14.70	38	Andrew Heiniger
12	Clock Ice Tea	16	25.20	Summary Count : 4	
13	Telephone Clock	23	54.00	Column Footer	
15	Ice Tea Chair	8	14.10	Page Footer Page 1 of 1	
18	Chair Clock	52	77.40		
24	Chair Shoe	21	21.60		
25	Shoe Shoe	23	24.30		
29	Chair Telephone	12	16.80		
36	Ice Tea Iron	16	4.80		
38	Clock Ice Tea	11	32.40		
40	Shoe Clock	33	8.40		
42	Iron Chair	13	29.70		
45	Iron Ice Tea	8	27.00		
48	Clock Clock	70	126.00		
Summary Count : 17					
Column Footer Total : 356 570.90					
Page Footer Page 1 of 1					

### New York

6 addresses

Products ordered by people in New York						Addresses in New York	
Title						Title	
Page Header						Page Header	
ID	Name	Qty	Price	ID	Name	Qty	Price
1	Chair Shoe	18	37.20	20	Ice Tea Telephone	7	30.00
2	Telephone Clock	15	37.20	21	Chair Chair	23	17.10
3	Chair Chair	18	38.10	22	Iron Iron	16	9.90
4	Ice Tea Shoe	56	96.00	23	Shoe Chair	2	11.40
5	Clock Clock	29	106.20	24	Chair Shoe	23	32.40
6	Ice Tea Chair	38	44.10	27	Telephone Iron	39	69.00
7	Telephone Shoe	49	37.80	28	Clock Iron	20	4.50
8	Ice Tea Clock	38	101.70	31	Ice Tea Telephone	13	7.20
10	Telephone Ice Tea	30	61.20	32	Clock Iron	26	57.00
11	Telephone Iron	23	26.40	33	Iron Chair	25	81.90
14	Telephone Iron	30	37.20	34	Chair Iron	17	76.80
15	Ice Tea Chair	50	56.40	35	Telephone Shoe	17	34.20
16	Ice Tea Shoe	27	58.20	38	Clock Ice Tea	31	97.20
17	Clock Ice Tea	23	33.00	40	Shoe Clock	52	12.60
18	Chair Clock	21	51.60	41	Clock Ice Tea	21	61.80
19	Ice Tea Ice Tea	37	33.00	43	Telephone Clock	29	28.20
Column Footer Total : 502 855.30						Summary Count : 6	
						Column Footer	
						Page Footer Page 1 of 1	
Column Footer Total : 863 1486.50							
Page Footer Page 1 of 2							

## City List

### New York (continued)

Page Header

ID	Name	Qty	Price
44	Ice Tea Ice Tea	58	72.00
45	Iron Ice Tea	37	81.00
46	Ice Tea Clock	12	9.30
47	Ice Tea Iron	10	26.70
48	Clock Clock	38	94.50
49	Iron Iron	31	6.60

Summary Count : 38

Column Footer Total : 1049 1776.60

Page Footer Page 2 of 2

### Olten 5 addresses

Title Products ordered by people in Olten

Page Header

ID	Name	Qty	Price	ID	Name	Qty	Price
0	Iron Iron	15	16.20	38	Clock Ice Tea	3	32.40
1	Chair Shoe	4	37.20	40	Shoe Clock	11	4.20
4	Ice Tea Shoe	13	19.20	41	Clock Ice Tea	41	61.80
5	Clock Clock	70	141.60	43	Telephone Clock	13	14.10
7	Telephone Shoe	13	25.20	45	Iron Ice Tea	25	54.00
8	Ice Tea Clock	22	33.90	46	Ice Tea Clock	22	18.60
9	Clock Telephone	18	25.80	47	Ice Tea Iron	26	53.40
10	Telephone Ice Tea	29	61.20	48	Clock Clock	22	31.50
11	Telephone Iron	24	13.20	49	Iron Iron	9	3.30
12	Clock Ice Tea	47	50.40				
13	Telephone Clock	8	27.00				
16	Ice Tea Shoe	8	29.10				
17	Clock Ice Tea	5	33.00				
18	Chair Clock	14	25.80				
19	Ice Tea Ice Tea	7	49.50				
20	Ice Tea Telephone	2	60.00				
21	Chair Chair	82	68.40				
22	Iron Iron	11	9.90				
23	Shoe Chair	20	11.40				
24	Chair Shoe	25	21.60				
25	Shoe Shoe	9	24.30				
28	Clock Iron	9	4.50				
30	Shoe Iron	14	34.80				
32	Clock Iron	21	28.50				
33	Iron Chair	17	27.30				
35	Telephone Shoe	15	11.40				
36	Ice Tea Iron	20	4.80				
37	Clock Shoe	24	14.10				

Column Footer Total : 566 909.30 Column Footer Total : 738 1182.60

Page Footer Page 1 of 2

Title Addresses in Olten

Page Header

ID	Name	Street
1	Susanne King	366 - 20th Ave.
24	James Sommer	333 Upland Pl.
30	Bill Sommer	362 - 20th Ave.
31	Bob Ringer	371 College Av.
49	George Fuller	534 - 20th Ave.

Summary Count : 5

Column Footer

Page Footer Page 1 of 1

## City List

### Olten (continued)

Summary Count : 37

### Oslo 6 addresses

Products ordered by people in Oslo							Addresses in Oslo			
Page Header							Page Header			
ID	Name	Qty	Price	ID	Name	Qty	Price	ID	Name	Street
0	Iron Iron	31	24.30	43	Telephone Clock	45	42.30	14	Bill Karsen	53 College Av.
1	Chair Shoe	24	111.60	44	Ice Tea Ice Tea	37	57.60	20	Janet Schneider	309 - 20th Ave.
2	Telephone Clock	35	74.40	45	Iron Ice Tea	32	81.00	26	James Clancy	195 Upland Pl.
3	Chair Chair	31	114.30	46	Ice Tea Clock	15	18.60	33	Mary King	491 College Av.
4	Ice Tea Shoe	41	57.60	47	Ice Tea Iron	33	80.10	42	Robert Ott	503 Seventh Av.
6	Ice Tea Chair	24	14.70					45	Janet May	396 Seventh Av.
7	Telephone Shoe	13	25.20					Summary		Count : 6
8	Ice Tea Clock	44	101.70					Column Footer		
9	Clock Telephone	23	51.60					Page Footer		Page 1 of1
10	Telephone Ice Tea	54	91.80							
11	Telephone Iron	31	39.60							
12	Clock Ice Tea	23	50.40							
13	Telephone Clock	17	27.00							
14	Telephone Iron	15	37.20							
15	Ice Tea Chair	24	28.20							
16	Ice Tea Shoe	18	58.20							
17	Clock Ice Tea	6	33.00							
18	Chair Clock	30	77.40							
19	Ice Tea Ice Tea	19	33.00							
20	Ice Tea Telephone	16	60.00							
21	Chair Chair	22	17.10							
23	Shoe Chair	45	45.60							
24	Chair Shoe	34	32.40							
25	Shoe Shoe	36	72.90							
26	Shoe Shoe	18	37.80							
27	Telephone Iron	18	69.00							
28	Clock Iron	36	9.00							
30	Shoe Iron	29	104.40							
31	Ice Tea Telephone	40	14.40							
32	Clock Iron	52	114.00							
33	Iron Chair	8	27.30							
35	Telephone Shoe	21	22.80							
36	Ice Tea Iron	20	4.80							
38	Clock Ice Tea	8	32.40							
39	Ice Tea Shoe	44	69.30							
40	Shoe Clock	38	21.00							
41	Clock Ice Tea	40	61.80							
42	Iron Chair	19	29.70							
Column Footer		Total :	1047 1896.90	Column Footer		Total :	1209 2176.50			
Page Footer							Page 1 of2			

## City List

### Oslo (continued)

Summary Count : 43

### Palo Alto 2 addresses

#### Title Addresses in Palo Alto

Page Header

ID	Name	Street
6	Laura White	506 Upland Pl.
12	George Ott	381 Upland Pl.

Summary Count : 2

Column Footer

Page Footer Page 1 of 1

### Paris 3 addresses

#### Title Products ordered by people in Paris

Page Header

ID	Name	Qty	Price
10	Telephone Ice Tea	12	30.60
20	Ice Tea Telephone	17	30.00
24	Chair Shoe	6	10.80
39	Ice Tea Shoe	16	23.10

Summary Count : 4

Column Footer Total : 51 94.50

Page Footer Page 1 of 1

#### Title Addresses in Paris

Page Header

ID	Name	Street
5	Laura Miller	294 Seventh Av.
18	Sylvia Fuller	158 - 20th Ave.
25	Sylvia Steel	269 College Av.

Summary Count : 3

Column Footer

Page Footer Page 1 of 1

### San Francisco 3 addresses

#### Title Addresses in San Francisco

Page Header

ID	Name	Street
3	Michael Clancy	542 Upland Pl.
7	James Peterson	231 Upland Pl.
48	Robert White	549 Seventh Av.

Summary Count : 3

Column Footer

Page Footer Page 1 of 1

### Seattle 5 addresses

#### Title Products ordered by people in Seattle

Page Header

ID	Name	Qty	Price	ID	Name	Qty	Price
0	Iron Iron	30	16.20	5	Clock Clock	18	35.40
1	Chair Shoe	60	148.80	7	Telephone Shoe	56	50.40
2	Telephone Clock	4	37.20	8	Ice Tea Clock	30	101.70
4	Ice Tea Shoe	16	19.20	9	Clock Telephone	38	77.40

Column Footer Total : 110 221.40 Column Footer Total : 252 486.30

Page Footer Page 1 of 2

#### Title Addresses in Seattle

Page Header

ID	Name	Street
8	Andrew Miller	288 - 20th Ave.
15	Bill Clancy	319 Upland Pl.
21	Julia Clancy	18 Seventh Av.
27	Bob Sommer	509 College Av.

Column Footer

Page Footer Page 1 of 2



## City List

### Seattle (continued)

Page Header

ID	Name	Qty	Price
10	Telephone Ice Tea	17	30.60
11	Telephone Iron	21	13.20
13	Telephone Clock	28	54.00
15	Ice Tea Chair	38	28.20
16	Ice Tea Shoe	3	29.10
17	Clock Ice Tea	45	66.00
18	Chair Clock	20	51.60
19	Ice Tea Ice Tea	25	33.00
20	Ice Tea Telephone	44	120.00
21	Chair Chair	8	17.10
23	Shoe Chair	18	11.40
24	Chair Shoe	12	10.80
26	Shoe Shoe	17	113.40
27	Telephone Iron	42	103.50
28	Clock Iron	18	4.50
29	Chair Telephone	17	16.80
30	Shoe Iron	44	139.20
31	Ice Tea Telephone	17	7.20
32	Clock Iron	14	28.50
33	Iron Chair	7	27.30
34	Chair Iron	16	38.40
35	Telephone Shoe	33	22.80
37	Clock Shoe	9	14.10
38	Clock Ice Tea	56	129.60
39	Ice Tea Shoe	17	23.10
40	Shoe Clock	16	8.40
41	Clock Ice Tea	24	61.80
42	Iron Chair	16	29.70
43	Telephone Clock	35	28.20
44	Ice Tea Ice Tea	19	14.40
46	Ice Tea Clock	40	27.90
49	Iron Iron	34	9.90

Summary

Count : 40

Column Footer Total : 1022 1800.00

Page Footer

Page 2 of 2

Page Header

ID	Name	Street
34	Julia May	33 Upland Pl.

Summary Count : 5

Column Footer

Page Footer

Page 2 of 2