
Address Report

ID	Name	Street
1. Berne		
22	Bill Ott	250 - 20th Ave.
9	James Schneider	277 Seventh Av.
		Count : 2
2. Boston		
32	Michael Ott	339 College Av.
23	Julia Heiniger	358 College Av.
		Count : 2
3. Chicago		
39	Mary Karsen	202 College Av.
35	George Karsen	412 College Av.
11	Julia White	412 Upland Pl.
		Count : 3
4. Dallas		
47	Janet Fuller	445 Upland Pl.
43	Susanne Smith	2 Upland Pl.
40	Susanne Miller	440 - 20th Ave.
36	John Steel	276 Upland Pl.
37	Michael Clancy	19 Seventh Av.
19	Susanne Heiniger	86 - 20th Ave.
10	Anne Fuller	135 Upland Pl.
4	Sylvia Ringer	365 College Av.
0	Laura Steel	429 Seventh Av.
		Count : 9
5. Lyon		
38	Andrew Heiniger	347 College Av.
28	Susanne White	74 - 20th Ave.
17	Laura Ott	443 Seventh Av.
2	Anne Miller	20 Upland Pl.
		Count : 4
6. New York		
46	Andrew May	172 Seventh Av.
44	Sylvia Ott	361 College Av.
41	Bill King	546 College Av.
29	Andrew Smith	254 College Av.
16	John Fuller	195 Seventh Av.
13	Laura Ringer	38 College Av.
		Count : 6

ID	Name	Street
7. Olten		
49	George Fuller	534 - 20th Ave.
30	Bill Sommer	362 - 20th Ave.
31	Bob Ringer	371 College Av.
24	James Sommer	333 Upland Pl.
1	Susanne King	366 - 20th Ave.
		Count : 5
8. Oslo		
45	Janet May	396 Seventh Av.
42	Robert Ott	503 Seventh Av.
33	Mary King	491 College Av.
26	James Clancy	195 Upland Pl.
20	Janet Schneider	309 - 20th Ave.
14	Bill Karsen	53 College Av.
		Count : 6
9. Palo Alto		
12	George Ott	381 Upland Pl.
6	Laura White	506 Upland Pl.
		Count : 2
10. Paris		
25	Sylvia Steel	269 College Av.
18	Sylvia Fuller	158 - 20th Ave.
5	Laura Miller	294 Seventh Av.
		Count : 3
11. San Francisco		
48	Robert White	549 Seventh Av.
7	James Peterson	231 Upland Pl.
3	Michael Clancy	542 Upland Pl.
		Count : 3
12. Seattle		
34	Julia May	33 Upland Pl.
27	Bob Sommer	509 College Av.
21	Julia Clancy	18 Seventh Av.
15	Bill Clancy	319 Upland Pl.
8	Andrew Miller	288 - 20th Ave.
		Count : 5

*There were 50 address records on this report from the following cities :
Berne, Boston, Chicago, Dallas, Lyon, New York, Olten, Oslo, Palo Alto, Paris, San Francisco,
Seattle.*

Hello! I'm the report's scriptlet object.