
Address Report

CustomBeanFactory.java - Bean Collection

ID	Name	Street
1. Berne		
9	James Schneider	277 Seventh Av.
22	Bill Ott	250 - 20th Ave.

Count : 2

2. Boston		
23	Julia Heiniger	358 College Av.
32	Michael Ott	339 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.

Count : 3

ID	Name	Street
7. Oslo		
45	Janet May	396 Seventh Av.
42	Robert Ott	503 Seventh Av.

Count : 2

8. Paris		
25	Sylvia Steel	269 College Av.
18	Sylvia Fuller	158 - 20th Ave.
5	Laura Miller	294 Seventh Av.

Count : 3

9. San Francisco		
48	Robert White	549 Seventh Av.
7	James Peterson	231 Upland Pl.

Count : 2

There were 30 address records on this report.