

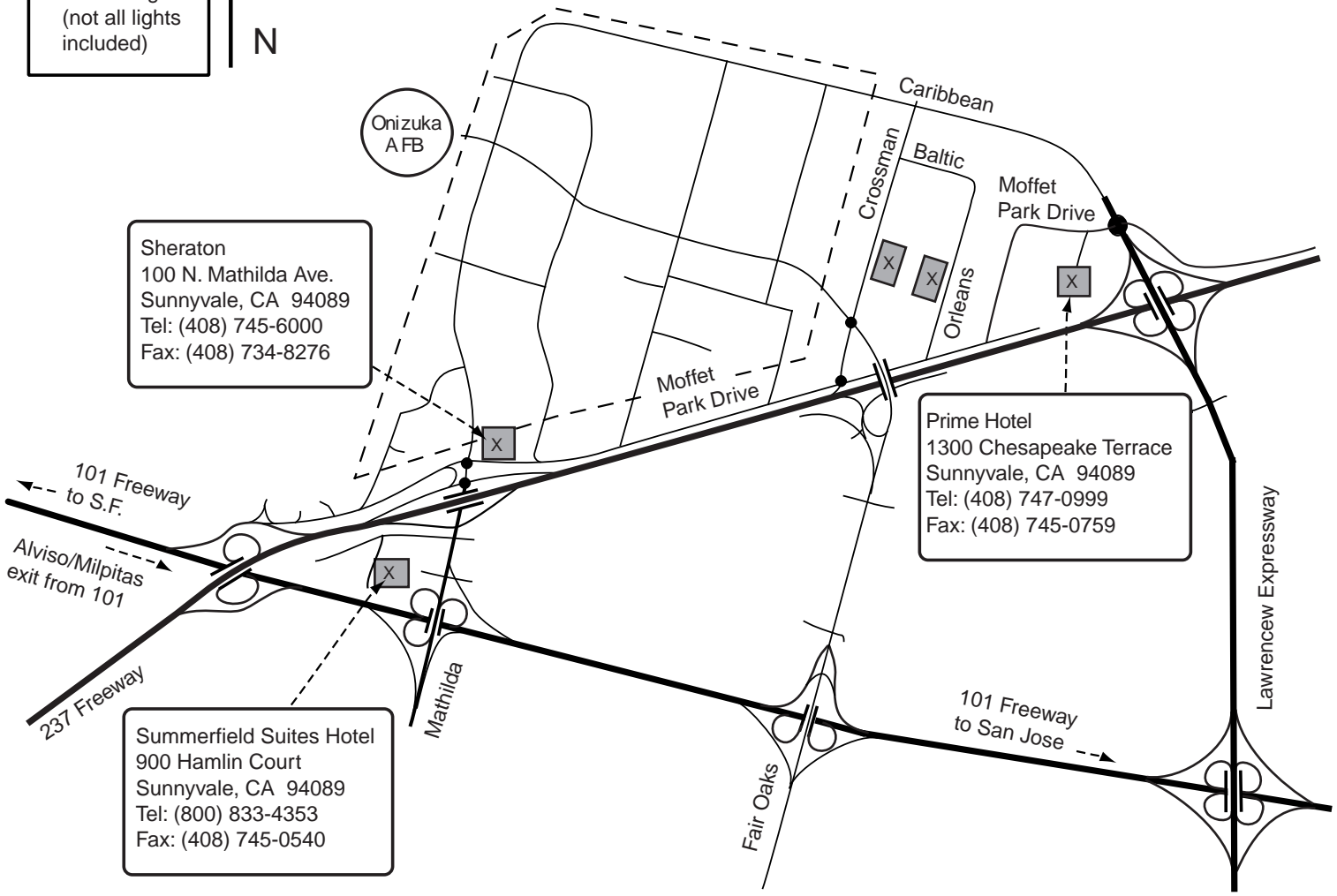
Map to Molecular Devices Corp

1312 Crossman Avenue
Sunnyvale, CA 94089
Tel: (408) 747-1700
Fax: (408) 747-3601

1311 Orleans Drive
Sunnyvale, CA 94089
Tel: (408) 747-1700
Fax: (408) 747-3601

● = traffic lights
(not all lights included)

↑ N



From San Francisco Airport:

To Crossman Building: Take US-101 South towards San Jose. Stay on 101 South for approximately 27 miles, then take Lawrence Expressway North. Continue on Lawrence Expressway until you cross over Highway 237 (at this point, Lawrence turns into Caribbean). Go past the first stop light after crossing Highway 237, then make a left onto Crossman Avenue. Molecular Devices is on the left side of the street at 1312 Crossman Avenue.

To Orleans Building: Take US-101 South towards San Jose. Stay on 101 South for approximately 27 miles, then take Lawrence Expressway North. Continue on Lawrence Expressway until you cross over Highway 237 (at this point, Lawrence turns into Caribbean). At the first stop light after crossing Highway 237, make a left onto Moffet Park Drive (note: the overhead sign will read "Baylands Park"). Follow Moffet Park Drive as it curves left, then right. Make a right onto Orleans Drive. Molecular Devices is on the left side of the street at 1311 Orleans Drive.

From San Jose Airport:

To Crossman Building: Make a left onto the Guadalupe Parkway (Highway 87). After approximately 3 miles, exit onto US-101 North. Take the Lawrence Expressway North exit. Continue on Lawrence Expressway until you cross over Highway 237 (at this point, Lawrence turns into Caribbean). Go past the first stop light after crossing Highway 237, then make a left onto Crossman Avenue. Molecular Devices is on the left side of the street at 1312 Crossman Avenue.

To Orleans Building: Make a left onto the Guadalupe Parkway (Highway 87). After approximately 3 miles, exit onto US-101 North. Take the Lawrence Expressway North exit. Continue on Lawrence Expressway until you cross over Highway 237 (at this point, Lawrence turns into Caribbean). At the first stop light after crossing Highway 237, make a left onto Moffet Park Drive (note: the overhead sign will read "Baylands Park"). Follow Moffet Park Drive as it curves left, then right. Make a right onto Orleans Drive. Molecular Devices is on the left side of the street at 1311 Orleans Drive.