

TRAVEL INFORMATION

HOTELS

This list is intended as a guide in locating a hotel near the training facility. The hotels below offer a corporate rate for employees and customers of Applied Biosystems. Please inquire with the hotel regarding the corporate rate.

Crown Plaza Hotel – Foster City (0.4 miles to AB)

Address: 1221 Chess Dr., Foster City, CA 94404

Phone: 650-570-5700

Fax: 650-570-0540

WWW: www.crowneplaza.com

Note: Complimentary shuttle to and from San Francisco International Airport.

Please inquire directly with the hotel regarding availability of a shuttle to and from Applied Biosystems.

Free high-speed internet in rooms.

Restaurants on site.

Hilton Garden Inn – San Mateo (0.7 miles to AB)

Address: 2000 Bridgepointe Cir., San Mateo, CA 94404

Phone: 650-522-9000

Fax: 650-522-9099

WWW: www.hiltongardeninn.com

Note: Complimentary shuttle to and from San Francisco International Airport.

Free high-speed internet in rooms.

Restaurant on site.

Hyatt Regency – San Francisco Airport (7.4 miles to AB)

Address: 1333 Bayshore Highway, Burlingame, CA 94010

Phone: 650-347-1234

Fax: 650-696-2669

WWW: www.sanfranciscoairport.hyatt.com

Note: Complimentary shuttle to and from San Francisco International Airport.

Parking fee \$18/day.

HSIA or Wireless \$9.95/day.

Restaurant on site.

Marriott Hotel – San Mateo (3.1 miles to AB)

Address: 1770 S. Amphlett Blvd., San Mateo, CA 94402

Phone: 650-653-6000

Fax: 650-653-6084

WWW: www.marriott.com

Note: Complimentary shuttle to and from San Francisco International Airport.
HSIA or Wireless \$9.95/day.

Restaurant on site.

Residence Inn – San Mateo (1.7 miles to AB)

Address: 2000 Winward Way, San Mateo, CA 94404

Phone: 650-574-4700

Fax: 650-572-9084

WWW: www.marriott.com/residenceinn

Note: Complimentary shuttle to and from Applied Biosystems.

Complimentary hot breakfast and light dinner.

Complimentary grocery shopping service.

Fully equipped kitchens.

Free high-speed internet in rooms.

Services provided by the hotels including shuttle information are subject to change at any time without notice, please confirm service availability with the hotel.

GROUND TRANSPORTATION

This information was obtained from the yellow pages and other public sources. Applied Biosystems is not affiliated with any company appearing on this list and accepts no responsibility for anyone who patronizes any of these businesses.

TAXI COMPANIES

Yellow Allied Cab

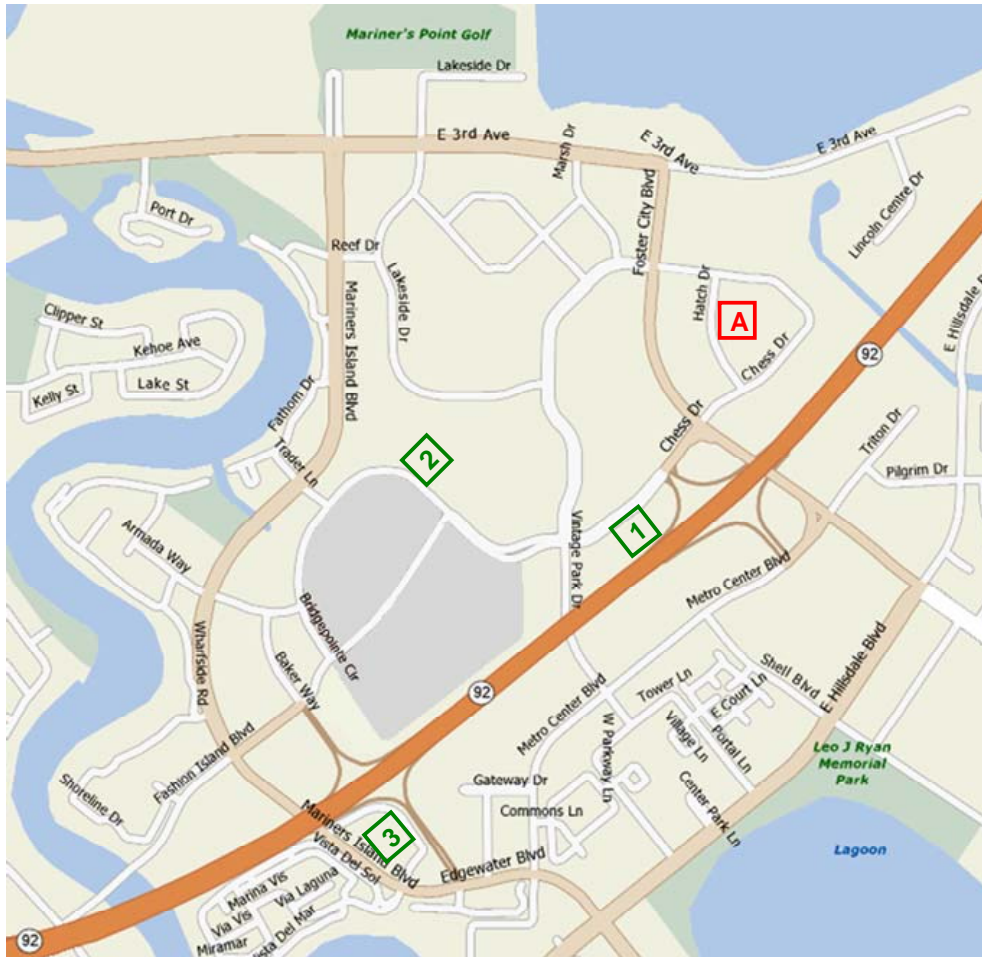
Phone: 650-854-4000
Toll-free: 800-220-1105
WWW: <http://www.yellowalliedcab.com>

Yellow Cab

Phone: 650-254-1234
Toll-free: 800-595-1222
WWW: <http://www.yellowcabpeninsula.com>

Yellow Cab AAA

Phone: 650-361-1234
Toll-free: 800-910-2227
WWW: <http://www.yellowcabaaa.com>



Applied Biosystems

A 353 Hatch Dr.
Foster City, CA 94404

Hotels

1 Crown Plaza Hotel
1221 Chess Dr.
650-570-5700

2 Hilton Garden Inn
2000 Bridgepointe Cir.
650-522-9000

3 Residence Inn
2000 Winward Way
650-574-4700