

DRIVING DIRECTIONS AND TRAVEL INFORMATION

From Logan Airport via Mass Pike to Applied Biosystems—Use the Ted Williams Tunnel, which will bring you directly to the westbound Massachusetts Turnpike (I-90). Proceed west on the Turnpike to Exit 13 (Natick).

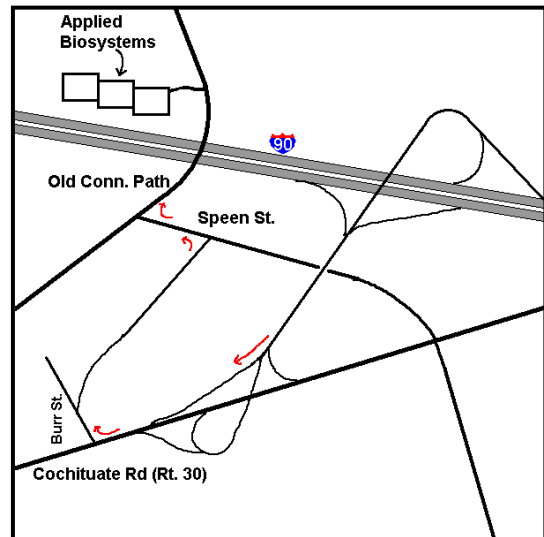
See map below right for remaining directions. Bear right as you leave the tollbooths and take the exit ramp onto Route 30 West. At the first light make a right onto Burr St. At Speen St. (light), make a left. At the next light make a right onto Old Connecticut Path (there is a right turn-off with stop sign). You will pass over the Massachusetts Turnpike, before coming to the Carousel Office Center on your left at 500 Old Connecticut Path.

From Route 128/95 North or South to Applied Biosystems—Take exit for Mass Pike (I-90) west. Proceed on the Massachusetts Turnpike to Exit 13 (Natick). Follow directions in the second paragraph above (starting “See map below right...”).

From Route 9 Westbound to Applied Biosystems—In Natick, take the exit for Mass Pike/Speen Street (If you get to the Natick Mall, you have gone too far). Stay in the right lane and take Speen St. across Route 30 (See map above right). Continue about ½ mile to the end of Speen Street and make a right at the light onto Old Connecticut Path. You will pass over the Mass Pike, before coming to the Carousel Office Center on your left at 500 Old Connecticut Path.

From Route 9 Eastbound to Applied Biosystems—Take the first exit to the right after the traffic lights at the Natick Mall. Almost immediately after exiting Route 9, there will be a secondary exit marked Speen Street – Mass Pike. (Be careful not to take the exit for Speen Street – Natick.) You will make a left over Route 9. As you cross over, get into the right lane and take Speen St. across Route 30 (See map above). Continue about ½ mile to the end of Speen Street and make a right at the light onto Old Connecticut Path. You will pass over the Mass Pike, before coming to the Carousel Office Center on your left at 500 Old Connecticut Path.

At 500 Old Connecticut Path—The main entrance to Applied Biosystems is located in Building B, the second building as indicated on the map. Visitors (not employees) may park in the visitor parking areas near the flagpoles; then proceed to the security desk.



TRAVEL INFORMATION

HOTELS

Hampton Inn

319 Speen St (Route 9 and Speen St.), Natick, MA 01760-1506

Phone: 508-653-5000

Fax: 508-651-9733

WWW: <http://www.hamptoninn.com>

Note: Complimentary continental breakfast. Government rates maybe available.

Crowne Plaza Boston-Natick

1360 Worcester Road (Route 9), Natick, MA 01760-1514

Phone: 508-653-8800

Fax: 508-653-170

<http://www.ihg.com/h/d/CP/1/en/rates?hotelCode=BOSNK&rateCode=ILP6P&IATAno=99502056>

Note: Complimentary shuttle service to and from Applied Biosystems and from the Logan Express Framingham terminal. Government rates maybe available.

Residence Inn Framingham

400 Staples Drive, Framingham, Massachusetts 01701

Phone: 508-370-0001

Fax: 508-370-0440

WWW: <http://marriott.com>

Note: A rental car is recommended if you plan to stay at this hotel.

Courtyard By Marriott

342 Speen St., Natick, MA 01760

Phone: 508-655-6100

Fax: 508-655-6104

Note: Local area shuttle upon request.

GROUND TRANSPORTATION

SHUTTLE (BUS) SERVICE

Logan Express Framingham

Phone: 800-235-6426

WWW: http://www.massport.com/logan/getti_typeo_logan.html

Fare: \$11.00 each way or \$20.00 round trip.

Note: This service will bring you to the Logan Express Framingham bus station.

The Crowne Plaza provides complimentary shuttle service from the bus station to the hotel (call the hotel upon arrival at the bus station).

TAXI COMPANIES

Natick Cab

Phone: 508-653-2300

Fare from Airport: \$49 one way; includes toll; tip not included

Credit: Cash or Major Credit Card

Tommy's Taxi

Phone: 508-875-3500

Fare from Logan: \$50 one way; does not include toll or tip

Credit: Cash or American Express Credit Card
