

# MassBioHub Site Visit Information



## Location

The MassBioHub Conference & Business Center is in the same building as the MIT Credit Union (look for the bank sign outside). Enter the lobby and take the elevator to the 5th floor. Our office is on the right. Please check in with our Office Manager, Stephanie, at reception and let her know who you're meeting from the Hub team.



## Technology Square Garage

**800 Technology Square  
Cambridge, MA 02139**

Parking is available in the Technology Square Garage, just steps from our building. Use the sliding doors at the southeast corner to return to your car. Scan the QR code on your Flash Parking ticket for access.



## Daily Garage Parking Rates

0 – 1 hour: \$26.00 • 1 – 2 hours: \$34.00  
2 – 3 hours: \$40.00 • 24 hours: \$54.00

## Accepted Forms of Payment

American Express, Discover, MasterCard, and Visa.  
MassBio does not validate parking.

## MassBioHub Staff



**Nate Gaul**  
*Operations Manager*  
(617) 675-9715  
Nate.Gaul@massbio.org



**Stephanie Haskell**  
*Office Manager*  
(617) 694-4193  
Stephanie.Haskell@massbio.org



**Emily Hill**  
*Sales Specialist*  
(617) 274-9633  
Emily.Hill@massbio.org



**Bailey Nascimento**  
*Events Manager*  
(617) 997-1675  
Bailey.Nascimento@massbio.org