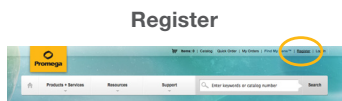




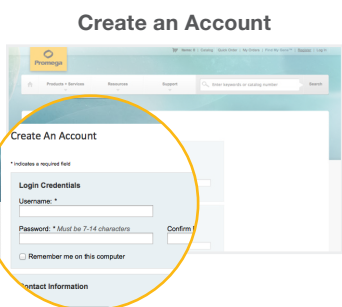
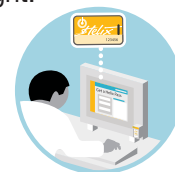
User Guide

Helix is the only on-site stocking program that automates the entire purchasing process—from obtaining products to invoicing, inventory and re-order—with the simple swipe of a card.



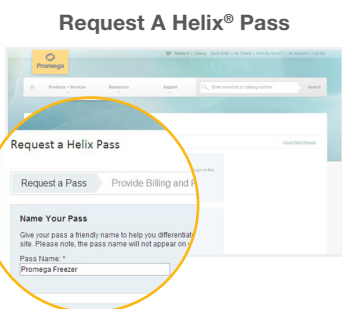
How to Register for Helix

- Visit www.promega.com and select “**Register**” on the top right.
- Follow the prompts to set up your promega.com profile.
- Click on any of the Helix® links on the account landing page.
- Follow the prompts to set up a pass.
- Download the mobile app from the iTunes App Store or Google Play.
- Your pass will arrive in 1–2 business days.



How to Access Helix

See what’s in your Helix® unit using the mobile app, touchscreen or promega.com.



Use the mobile app, touchscreen or pass to unlock Helix.



- Remove your product and close the door.
- You will receive an order confirmation e-mail.
- Log into your account at www.promega.com to review your order history.

Problems or questions?

Please contact the Promega Helix® team at helix@promega.com or 1-800-356-9526, Extension 2222 for assistance.

If you need a product that is not in your unit, e-mail your request to helix@promega.com

