



The products you need. . . Exactly when you need them!

To access PromegaExpress, you need a pass. Sign up on-line:

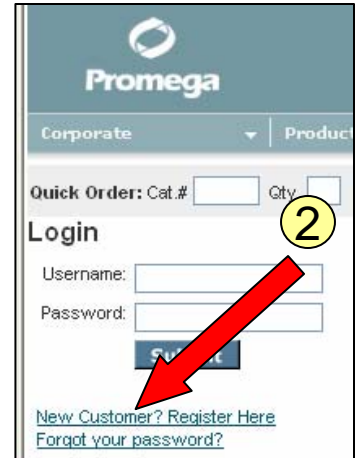
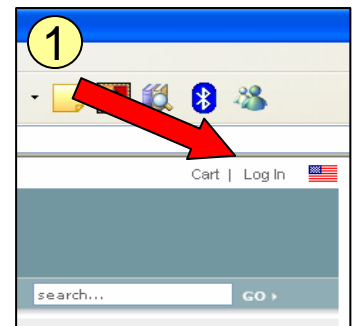
You will need:

1. Your PromegaExpress site-code
2. A valid purchase order number or credit card
3. Your contact information, including mailing address, so we can send you your pass



How to register for a pass on-line:

1. Go to promega.com and select “Login” from the upper right corner by your country’s flag icon.
2. On the left side of the screen under the Login fields click on “[New customer? Register here](#)”.
3. Complete the form with your information. You must have your site-code to receive a pass, this can be obtained using the site-locator tool or from your Promega Representative. You may also need the "Promega Partnership Program" (PPP) code (not all freezers have a PPP code). If you need a PPP code it will be provided with your site-code. Click "Save".



PromegaExpress™ Program
 PromegaExpress™ Site Code:
[Promega Express™ Program Information](#)

Promega Partnership Program:
 Promega Partnership Program:
[Promega Partnership Program Information](#)

3

4. Enter your preferred billing method and valid purchase order or credit card information and click “Save”.
5. You will receive a confirmation email within a few minutes from USExpress@promega.com, **Keep this email**. Your pass will arrive in 1-2 days. To activate your pass reply to the email message.

Credit Card

Name on card:
 Card number:
 Type:
 Expiration month:
 Expiration year:

Purchase Order

Purchase Order Number:
 Saved PO Numbers:

Internal Reference Number:
 Tax information (if applicable to this account)
 Yes, please apply tax exempt status to this order. If we do not currently have your tax-exempt certificate on file, please fax it to 1-800-356-1970.
 Tax Exempt Certificate #:

4