

Product Number  
**826-05213**

## KB<sup>Plus</sup>™ 10X TBE Buffer Pack

### Description

KB<sup>Plus</sup> 10X TBE buffer pack contains enough pre-measured powder to make one liter of 10X TBE. The buffer is optimized for use with KB<sup>Plus</sup> Gel Matrix products. The 10X stock solution contains 0.89 M Tris, 0.89 M borate, and 0.02 M EDTA. KB<sup>Plus</sup> products are specifically designed to optimize read length, accuracy, and run time with LI-COR automated infrared DNA analyzers.

Quantity: 1 pack

Storage: Ambient

Revised: September 2012

Updates available at:  
<http://biosupport.licor.com>

### Protocol

#### To prepare 10X TBE:

Pour the contents of the 10X TBE Buffer Pack into a one-liter beaker. Add enough distilled water (18 MΩ) to bring volume to about 800 mL. Stir the solution well until all of the solids have gone into solution and the solution is clear. Transfer beaker contents to a one-liter graduated cylinder, and add distilled water to bring the final volume to 1000 mL. Store the 10X TBE stock solution at room temperature, and use within three months.

#### To prepare 0.8X gel running buffer:

Add 80 mL of 10X TBE to 920 mL of distilled water (18 MΩ). Mix well.

#### To prepare 1.0X gel running buffer:

Add 100 mL of 10X TBE to 900 mL of distilled water (18 MΩ). Mix well.

### Storage

Store dry buffer pack at room temperature. Expiration is one year from date received.

© 2012 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR and KB<sup>Plus</sup> are trademarks and registered trademarks of LI-COR, Inc. in the United States and other countries.

Doc #988-13238



4647 Superior Street • P.O. Box 4000  
Lincoln, Nebraska 68504 USA  
North America: 800-645-4267  
International: 402-467-0700  
FAX: 402-467-0819

LI-COR GmbH Germany, Serving Europe,  
Middle East and Africa: +49 (0) 6172 17 17 771  
LI-COR UK Ltd. UK, Serving UK, Ireland, and  
Scandinavia: +44 (0) 1223 422104  
All other countries, contact LI-COR Biosciences  
or a local LI-COR distributor:

<http://www.licor.com/distributors>

[www.licor.com/bio](http://www.licor.com/bio)

