

Product Number  
**827-05214**

## KB<sup>Plus</sup>™ 3.7% Gel Matrix

### Description

KB<sup>Plus</sup> Gel Matrix products are specifically designed to optimize read length, accuracy, and run time with LI-COR automated infrared DNA analyzers. The 3.7% Gel Matrix is designed for sequencing with 66 cm, 0.2 mm gel apparatus. Each bottle contains 3.7% KB<sup>Plus</sup> modified acrylamide, 7 M Urea, and 1X TBE buffer.

Quantity: 500 mL

Storage: 4 °C

Revised: September 2012

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

### Protocol

#### To prepare 3.7% gel solution for use with 66 cm, 0.2 mm gels:

Bring 40 mL KB<sup>Plus</sup> 3.7% Gel Matrix to room temperature (10-15 minutes) in a beaker. During that time, prepare glass plates for gel injection. When ready to inject gel solution, add 230 µL 10% APS, and 23 µL TEMED. Mix to homogenate and inject into gel cassette. After injection, allow at least 1.5 hour for polymerization before use.

Pre-run gel for 30 minutes (0.8X TBE running buffer) using the following electrophoresis conditions:

Voltage	3000 V
Current	30 mA
Power	80 W
Temperature	45 °C

After the pre-run, load samples and begin electrophoresis.

Expiration: Six months from receipt.

© 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 #988-13239



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)