

Product Number
827-05215

KB^{Plus}™ 5.5% Gel Matrix

Description

KB^{Plus} Gel Matrix products are specifically designed to optimize read length, accuracy, and run time with LI-COR automated infrared DNA analyzers. The 5.5% Gel Matrix is designed for sequencing with 41 cm, 0.2 mm gel apparatus. Each bottle contains 5.5% KB^{Plus} 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 5.5% gel solution for use with 41 cm, 0.2 mm gels:

Bring 30 mL KB^{Plus} 5.5% 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 200 µL 10% APS and 20 µL TEMED. Swirl gently to homogenate and inject into gel cassette. After injection, allow at least 1.5 hour for polymerization before use.

Pre-run gel for 15 minutes (0.8X TBE running buffer) using the following electrophoresis conditions:

Voltage	2000 V
Current	25 mA
Power	50 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^{Plus} are trademarks and registered trademarks of LI-COR, Inc. in the United States and other countries.

Doc #988-13240



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

