

Product Number
827-05607

KB^{Plus}™ 6.5% Gel Matrix for Genotyping

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 6.5% Gel Matrix is designed for genotyping or AFLP[®] technology (Amplified Fragment Length Polymorphism) with 25 cm, 0.25 mm gel apparatus. Each bottle contains 6.5% KB^{Plus} modified acrylamide, 7M Urea, and 1X TBE buffer.

Quantity: 500 mL

Storage: 4 °C

Revised: September 2012

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

Protocol

To prepare 6.5% gel solution:

1. Determine the thickness of gel to be poured (0.25 or 0.4 mm).
2. Using the table below, bring the correct volume of KB^{Plus} 6.5% Gel Matrix to room temperature (10-15 minutes). During that time, prepare glass plates for gel injection. When ready to inject gel solution, add the amounts of APS and TEMED indicated in the table. Swirl gently to homogenate and inject into gel cassette. After injection, center and insert the gel casting comb and allow at least 1.5 hour for polymerization before use.
3. Prerun gel for 10 minutes (KB^{Plus} 1X TBE running buffer, LI-COR P/N 826-05606) using the following electrophoresis conditions:
Voltage 1500 V
Current 40 mA
Power 40 W
Temperature 50 °C
4. After the prerun, load samples and begin electrophoresis.

Gel Thickness	KB ^{Plus}	APS	TEMED
0.25 mm	20 mL	150 µL	15 µL
0.40 mm	40 mL	300 µL	30 µL

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. AFLP is a registered trademark of KeyGene N.V.

Doc #988-13241



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