

Product Number
928-40012

Quantity: 2-pack

Storage: Ambient

Revised: September 2012

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

Tris Glycine Dry Powder Buffer

Description

A single strength (1X) working solution contains 0.025 M Tris base and 0.192 M glycine.

Appearance

Uniform white powder.

Protocol

Dissolve the contents of one pack in 1 liter of deionized water to make a 10X stock solution.

Storage

Store dry buffer packs in a tightly sealed container, at room temperature. Protect from moisture. Expiration is one year from date received.

© 2012 LI-COR, Inc. LI-COR is an ISO 9001 registered company. LI-COR is registered trademark of LI-COR, Inc. in the United States and other countries.

Doc #988-13206



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