

Novex™ WedgeWell™ Tris-Glycine Gels

Pub. No. MAN0014610 Rev. A.0

Instructions for performing electrophoresis using the Mini Gel Tank or the XCell SureLock™ Mini-Cell to are described below. For videos and detailed instructions visit **thermofisher.com**

Prepare Samples

Reagent	Reduced	Non-reduced
Sample	x μ L	x μ L
Tris-Glycine SDS Sample Buffer (2X)	20 μ L	20 μ L
NuPAGE™ Reducing Agent (10X)	4 μ L	—
<u>Deionized Water</u>	<u>to total volume</u>	<u>to total volume</u>
Total Volume	40 μ L	40 μ L

Heat samples at 85°C for 2 minutes.

Prepare 1X Buffer

Mini Gel Tank: For each chamber, mix 40 mL of 10X Novex™ Tris-Glycine SDS Running Buffer with 360 mL of deionized water.

XCell SureLock™ Mini-Cell: Mix 100 mL of 10X Novex™ Tris-Glycine SDS Running Buffer with 900 mL deionized water.

Load Sample

Load the appropriate concentration of your protein sample on the gel.

Run	Mini Gel Tank	
Conditions	Voltage	225V
	Run Time**:	25-40 minutes
	Expected Current*:	85–125 mA (start); 30–55 mA (end)

XCell SureLock™ Mini-Cell

Voltage	225V
Run Time**:	35-45 minutes
Expected Current*:	45–60 mA (start); 30–45 mA (end)

* Double the current setting for running two mini-gels.

** Run times may vary depending upon your power supply.

Important licensing information:

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/terms.and.conditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

© 2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit thermofisher.com/support

15 September 2015

ThermoFisher
S C I E N T I F I C