



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

MES SDS Buffer Kit (for Bis-Tris Gels)

For SDS Polyacrylamide Gel Electrophoresis
(Cat. #786-554)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

MES SDS Buffer Kit is a complete buffer kit to resolve small- to medium-sized proteins under denaturing conditions on Bis-Tris gels.

ITEM(S) SUPPLIED

Description	Cat. #786-554
MES SDS Running Buffer [20X]	2 x 250 ml
DTT [1M](15.4mg)	1 vial
Protein Antioxidant [200X]	15 ml
LDS Sample Loading Buffer [4X]	10 ml

STORAGE CONDITIONS

The kit is shipped at ambient temperature. Store at 4°C.

PREPARATION BEFORE USE

1. Add 100 µl deionized water to DTT vial. Mix well to dissolve and store it at -20°C after use.
2. Add 50 ml of MES SDS Running Buffer [20X] to 950 ml of deionized water to obtain 1X MES SDS Running Buffer.

PROTOCOL

Loading Sample Preparation

1. Bring the LDS Sample Loading Buffer [4X] to room temperature before use.
2. Prepare reduced or non-reduced protein samples as below:

Reagent	Reduced Protein Sample	Non-reduced Protein Sample
Protein Sample	x µl	x µl
LDS Sample Loading Buffer [4X]	2.5 µl	2.5 µl
DTT [0.5 M,10X]	0.5 µl	-

Make up the final volume of sample to 10 µl with deionized water.

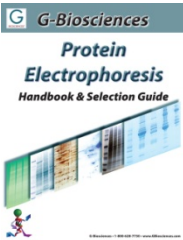
3. Vortex the tube to mix the contents.
4. Heat both the reduced and non-reduced samples at 70°C for 10 minutes and let it cool.
5. Centrifuge the sample tubes and load the samples on gel to run SDS-PAGE.

SDS-PAGE with Bis-Tris Gels and MES SDS Running Buffer

1. For reducing SDS-PAGE add 1 ml Protein Antioxidant [200X] to 200 ml of 1X MES SDS Running Buffer and fill the upper chamber of electrophoresis unit.
Note: *Reducing and non-reducing samples are preferably run in different gels. For non-reducing sample run, Protein Antioxidant is not added to the running buffer in the cathodic chamber of electrophoresis unit. If reducing and non-reducing samples are run on same gel for some reason, then do not use Protein Antioxidant.*
2. Load the samples and perform electrophoresis at constant voltage of 200 V (Approx. run time for gel is 35 min).

RELATED PRODUCTS

Download our Protein Electrophoresis Handbook.



<http://info2.gbiosciences.com/complete-protein-electrophoresis-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 1/30/2017 CMH



www.GBiosciences.com