



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Tris-Acetate SDS Buffer Kit (for Tris-Acetate Gels)

For SDS Polyacrylamide Gel Electrophoresis

(Cat. #786-589)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Tris-Acetate SDS Buffer Kit is a complete buffer kit to resolve high molecular weight proteins (36-400 kDa) under denaturing conditions on Tris/Acetate gels

## ITEM(S) SUPPLIED

Description	Cat. #786-589
Tris-Acetate/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 Tris-Acetate/SDS Running Buffer to 950 ml of deionized water to obtain 1X Tris-Acetate/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 [ 1M](15.4mg)	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 Tris-Acetate Gels and Tris-Acetate/SDS Running Buffer

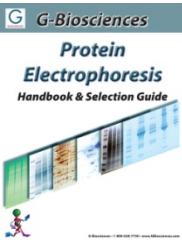
1. For reducing SDS-PAGE add 1 ml Protein Antioxidant [200X] to 200 ml of 1X Tris-Acetate/SDS Running Buffer and fill the upper chamber of electrophoresis unit.

**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 150 V (Approx. run time for gel is 1 hr).

### 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](http://www.GBiosciences.com) or contact us.

Last saved: 1/30/2017 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)