



431PR-02

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

A Geno Technology, Inc. (USA) brand name

SDS-PAGE Sample Loading Buffer

2X & 6X Concentrated Buffers For Running SDS
Polyacrylamide Gel Electrophoresis

(Cat. # 786-025 & 786-701)



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

INTRODUCTION

The SDS-PAGE Sample Loading Buffers are suitable for loading protein samples on to the SDS-polyacrylamide gels. The buffers are provided in 2X and 6X concentrations containing Tris-HCl, glycerol, SDS and bromophenol blue (BPB) in recommended concentrations and is stable at room temperature.

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-025	SDS-PAGE Sample Loading Buffer [2X]	2.5ml
786-701	SDS-PAGE Sample Loading Buffer [6X]	25ml

STORAGE CONDITIONS

Shipped at ambient temperature. Store at room temperature upon arrival.

ADDITIONAL ITEMS REQUIRED

- Reducing agent: β -mercaptoethanol or DTT

PREPARATION BEFORE USE

The SDS may precipitate at cooler temperatures. Simply warm in a 37°C waterbath or incubator to redissolve the SDS.

2X SDS-PAGE SAMPLE LOADING BUFFER PROTOCOL

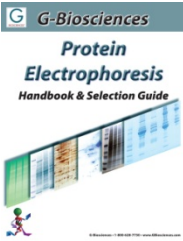
1. Add an equal volume of SDS-PAGE Sample Loading Buffer [2X] to the tube containing protein solution.
2. For reducing gels, add reducing agent to a final concentration of 2-5% β -mercaptoethanol or 5-20mM DTT.
3. Vortex the tube to mix the contents. Place the sample tube in a boiling water bath for 5-10 minutes.
4. After the boiling is complete, vortex and centrifuge the tube for 30 seconds. The sample is now ready for loading on SDS-PAGE gels. Vortex the tube before loading the protein solution on the gel.

6X SDS-PAGE SAMPLE LOADING BUFFER PROTOCOL

1. Add one volume of SDS-PAGE Sample Loading Buffer [6X] to five volumes of protein solution. For example add 5µl SDS-PAGE Sample Loading Buffer [6X] to 25µl protein solution.
2. For reducing gels, add reducing agent to a final concentration of 2-5% β-mercaptoethanol or 5-20mM DTT.
3. Vortex the tube to mix the contents. Place the sample tube in a boiling water bath for 5-10 minutes.
4. After the boiling is complete, vortex and centrifuge the tube for 30 seconds. The sample is now ready for loading on SDS-PAGE gels. Vortex the tube before loading the protein solution on the gel.

RELATED PRODUCTS

Download our Protein Electrophoresis Handbook.



<http://info.gbiosciences.com/complete-protein-sample-preparation-handbook/>

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

Last saved: 3/27/2014 CMH



www.GBiosciences.com