



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

A Geno Technology, Inc. (USA) brand name

LDS Sample Loading Buffer [4X]

For SDS Polyacrylamide Gel Electrophoresis

(Cat. #786-323)



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

ITEM(S) SUPPLIED

| Cat. # | Description | Size |
|---------|--------------------------------|-------|
| 786-323 | LDS Sample Loading Buffer [4X] | 10 ml |

STORAGE CONDITIONS

Shipped at ambient temperature. Store at 4°C.

ADDITIONAL ITEMS REQUIRED

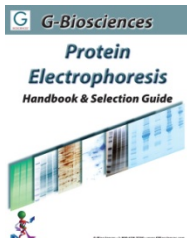
- Reducing agent: β -mercaptoethanol (Cat. # BC98), DTT (Cat. # RC-046) or TCEP.HCl (Cat. # 786-873)

PROTOCOL

- Bring the LDS Sample Loading Buffer [4X] to room temperature before use.
- Add one volume of LDS Sample Loading Buffer [4X] to three volumes of protein solution. For example, add 5 μ l LDS Sample Loading Buffer [4X] to 15 μ l protein solution.
- For reducing gels, add reducing agent to a final concentration of 2-5 % β -mercaptoethanol or 50 mM DTT.
- Vortex the tube to mix the contents.
- Heat the samples at 70°C for 10 minutes and let cool.
- Centrifuge the sample tubes and load the samples on gel to run SDS-PAGE.

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: 8/29/2016 CMH

This page is intentionally left blank



www.GBiosciences.com