



429PR-01

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

A Geno Technology, Inc. (USA) brand name

# SDS Detection & Estimation

(Cat. # 786-129)



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

## INTRODUCTION

SDS Detection & Estimation Kit is provided with reagents for the detection and estimation of SDS (sodium dodecyl sulfate) in a sample containing proteins or other agents. The working range of the assay is 0.02-0.2% SDS (1-10 $\mu$ g SDS).

## ITEM(S) SUPPLIED (Cat. # 786-129)

Description	Size
Blue Dye	30ml
Dye Extraction Buffer	15ml

## STORAGE CONDITIONS

The kit is shipped at ambient temperature. Upon arrival, store at room temperature.

## ADDITIONAL ITEMS NEEDED

- Chloroform,
- 15ml centrifuge tubes
- 1% SDS solution for standards

## PROTOCOL: SDS DETECTION

1. For SDS estimation, prepare an appropriate calibration plot with known concentrations of SDS ranging from 0.02% to 0.2%.

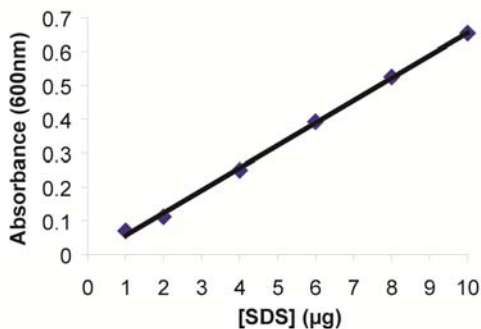


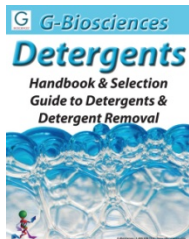
Figure 1: SDS Standard Curve

2. In a 15ml centrifuge tube, add 2ml Blue Dye and 1ml Dye Extraction Buffer for each sample.
3. Aliquot 1-5 $\mu$ l test sample or 5 $\mu$ l deionized water (blank) or 5 $\mu$ l Standards to the 15ml centrifuge tube. Vortex for 30 seconds.
4. Add 2ml chloroform into the mix, and mix the tube content by vortexing for 30 seconds.

5. Allow the tube to stand at room temperature for 5 minutes. If the lower, chloroform layer extracts color then it is indicative of the presence of SDS in the test sample.
6. Blank the spectrophotometer and then measure the optical density of the chloroform layer at 600nm. Ensure to use cuvettes that are compatible with chloroform, for example quartz cuvettes.

## RELATED PRODUCTS

Download our Detergents Handbook



<http://info.gbiosciences.com/complete-detergent-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.

Last saved: 7/26/2012 CMH



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