



SDS Detection & Estimation Kit

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 sensitivity is up to 2ng SDS present.

ITEM(S) SUPPLIED

Cat # 786-129

<i>DetergentOUT™</i> Blue Dye	30ml
<i>DetergentOUT™</i> Dye Extraction Buffer	15ml

STORAGE CONDITIONS

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

ITEMS NEEDED BUT NOT SUPPLIED

Chloroform, Centrifuge & Collection Tubes

PROTOCOL

DETECTION & ESTIMATION OF SDS

Mix 1-5µl of test sample with 2ml Blue Dye and 1ml Dye Extraction Buffer.

Prepare a control containing 1-5µl of water. [For estimation of SDS concentration, prepare a calibration plot using known concentrations of SDS]

Add 2ml chloroform into the mix, and mix the tube content by inverting the tube 5-6 times.

Allow the tube to stand at room temperature for 5 minutes. If the chloroform layer extracts color then it is indicative of the presence of SDS in the test sample. Measure the optical density of the chloroform layer at 600nm.

Compare the optical density of the chloroform layers of the control and the test sample.

Estimation of SDS Concentration:

Prepare an appropriate calibration plot with known concentrations of SDS. Use the calibration plots for the estimation of SDS concentration.

RELATED PRODUCTS

1. *Spin-OUT™*: Spin columns that are suitable for buffer exchange or removal of small molecules from protein and nucleic acid solutions.
2. *PROTEIN-Concentrate™ Kit*: concentrates dilute protein solution in a few minutes.
3. *Non-Interfering Protein Assay™*: A protein assay that is not affected by the presence of common laboratory agents such as detergents, reducing agents, EDTA, dyes etc.
4. *Tube-O-DIALYZER™*: Patented dialysis tube for dialyzing small samples.

NOTE: For other related products, visit our web site at www.GBiosciences.com or contact us.

Rev10.12.09/IA

