





# SDS PAGE Sample Buffer - 2x conc.

Cat. No.	Amount
BU-126	3 x 10 ml

### For general laboratory use.

Shipping: shipped at ambient temperature Storage Conditions: store at 4 °C Shelf Life: 12 months Form: liquid Concentration: 2x

### Description:

SDS PAGE Sample Buffer is the most commonly used sample buffer for Sodium Dodecyl Sulfate - Polyacrylamide Gel Electrophoresis (SDS-PAGE) of denatured proteins. SDS PAGE Sample Buffer ensures optimal band resolution when preparing proteins for SDS-PAGE with Tris-glycine-SDS running buffer.

### Content:

2x concentration of SDS PAGE Sample Buffer contains: 0.15 M Tris-HCl pH 6.8 10 % Mercaptoethanol 1.2 % SDS 30 % Glycerol 0.04 % Bromphenol blue

## Quality:

Molecular biology grade

### This mixture contains dangerous substances



Signal word: Danger

#### Hazard statements

H302 + H312 + H332 Harmful if swallowed, in contact with skin and if inhaled. H410 Very toxic to aquatic life with long lasting effects.

#### **Precautionary statements**

P261 Avoid breathing dust/fume/gas/mist/vapours/spray. P273 Avoid release to the environment. P280 Wear protective gloves/protective clothing/eve protection/fa

P280 Wear protective gloves/protective clothing/eye protection/face protection.

P301 + P312 IF SWALLOWED: Call a POISON CENTER/doctor/./ if you feel unwell.

P302 + P352 IF ON SKIN: Wash with plenty of soap and water. P304 + P340 IF INHALED: Remove person to fresh air and keep comfortable for breathing.

For further information see Material Safety Data Sheet.

