

9620 Medical Center Drive, Suite 101 Rockville, MD 20850, USA Web: www.abpbio.com

Product Information

5×SDS-PAGE Sample Loading Buffer

Catalog Number	Packaging Size
P013	5 mL

Storage upon receipt: • -20°C

Product Description

SDS-PAGE Sample Loading Buffer is a 5X solution of 250 mM Tris-HCl, pH 6.8, 10% SDS, 30% (v/v) Glycerol, 10 mM DTT, 0.05% (w/v) Bromophenol Blue for use in SDS-polyacrylamide gel electrophoresis of proteins.

Application

1. Warm the **SDS-PAGE Sample Loading Buffer (5X)** to room temperature.

2. Mix 4 volumes of protein sample with 1 volume of **SDS-PAGE Sample Loading Buffer (5X)**. Vortex briefly.

- 3. Heat to 95°C for 5 min to denature protein.
- 4. Cool to room temperature, and load on the gel.
- 5. Run electrophoresis until the blue dye reaches the bottom of the gel.