

Version: 02

Update: 09/30/2022

**Product name: 4X LDS Sample Buffer**

**Cat. No.: M00676-10/M00676-250**

**Description**

4X LDS Sample Buffer is used to prepare protein samples for denaturing polyacrylamide gel electrophoresis (PAGE) with GenScript's PAGE Gels and most other types of Bis-Tris gels. LDS sample buffer contains lithiumdodecyl sulfate with pH at 8.4, which helps reducing the disulfide bonds and ensure optimal protein separation.

**Note:** 4X LDS Sample Buffer is a highly viscous solution due to high concentration of LDS and glycerol. The buffer should be brought to room temperature (25°C) and mix well prior to use.

**General Specifications**

---

Buffer Type:	4X LDS Sample Buffer
Form:	Liquid
Product Size:	10 ml/250ml
Gel Compatibility:	GenScript's PAGE Gels, Most Bis-Tris PAGE Gels
Concentration(1X):	141 mM Tris Base, 106 mM Tris HCl, 2% LDS, 10% Glycerol, 0.51mM EDTA 0.22 mM SERVA Blue G-250, 0.175mM Phenol Red

---

## Preparing Samples

Follow the Instructions below to prepare samples for denaturing gel.

Reagent	Reduced Sample	Non-reduced Sample
Sample	x $\mu$ l	x $\mu$ l
4X LDS Sample Buffer	2.5 $\mu$ l	2.5 $\mu$ l
Reducing Agent 10x (500mM DTT or 25% $\beta$ -ME)	1 $\mu$ l	-
Deionized Water	Up to 6.5 $\mu$ l	Up to 7.5 $\mu$ l
Total Volume	10 $\mu$ l	10 $\mu$ l

Either  $\beta$ -ME or DTT can be used in the 4X LDS Sample Buffer. Make sure that a fresh solution of  $\beta$ -ME or DTT is used.

Heat samples at 70°C for 10 minutes, spin down the sample before loading.

**For best results, please add the reducing agent immediately prior to electrophoresis.**

## Storage

4°C for 12 months

**For research and manufacturing use. Direct human use, including taking orally and injection are forbidden.**

生产商：南京金斯瑞生物科技有限公司 江苏省南京市江宁区科学园雍熙路 28 号

Manufacturer: Nanjing GenScript Biotech Co., Ltd. No. 28 Yongxi Road, Jiangning District, Nanjing, Jiangsu, China