

Version: 02

Update: 09/30/2022

Product name: 5X Sample Buffer

Cat No: MB01015

Description

5X Sample Buffer is used as a tracking dye for SDS PAGE gel loading. This buffer contains SDS and is suitable for denaturing gel electrophoresis. Dilute 1:4 with sample before loading.

Note: 5X Sample Buffer should be brought to room temperature prior to use. Mix well before loading gel.

General Specifications

Buffer Type:	Sample Loading Buffer
Form:	Buffer
Concentrated:	5X
Product Size:	5 ml
Gel Compatibility:	PAGE Gels
Regulatory Statement:	Not for use in diagnostic procedures.

Preparing Samples

Follow the Instructions below to prepare samples for denaturing gel.

Reagent	Volume
Protein sample	x μl
5 x Sample Buffer	2µl
Deionized Water	Up to 8µl
Total Volume	10µl

Heat samples at 100°C for 10 minutes, spin down the sample before loading.

Contents and Storage

Contents: 5X Sample Buffer **Storage:** -20°C for 24 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