

Product Information

Catalog #	Description
Premixed Sample Buffers	
161-0737	Laemmli Sample Buffer , 30 ml
161-0738	Native Sample Buffer , 30 ml
161-0739	Tricine Sample Buffer , 30 ml
161-0767	Nucleic Acid Sample Buffer, 5x , 10 ml
161-0768	TBE-Urea Sample Buffer , 30 ml
161-0763	IEF Sample Buffer , 30 ml
161-0764	Zymogram Sample Buffer , 30 ml

Premixed Buffers

161-0744	10x Tris/Tricine/SDS , 1 L
161-0760	10x Tris/Tricine/SDS , 6 x 1 L
161-0734	10x Tris/Glycine , 1 L
161-0771	10x Tris/Glycine , 5 L

Reducing Agents

161-0710	2-mercaptoethanol , 25 ml
161-0610	Dithiothreitol , 1 g
161-0611	Dithiothreitol , 5 g

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Dr., Hercules, CA 94547 USA

510-741-1000

4006030 Rev E

Tricine Sample Buffer

Catalog #

161-0739

BIO-RAD

Introduction

Bio-Rad's tricine sample buffer is based on the method of Schagger von Jagow (1987) with modifications specially formulated in our laboratory to improve band tightness. The use of tricine sample buffer ensures optimal band resolution when preparing peptides and small proteins for SDS-PAGE with Tris-Tricine-SDS running buffer.

Specifications

Composition	200 mM Tris-HCl, pH 6.8 40% glycerol 2% SDS 0.04% Coomassie Blue G-250
Storage	Ambient temperature
Shelf life	1 year

Instructions for Use

1. Add Reducing Agent

Add 20 μl of β -mercaptoethanol per 950 μl of sample buffer for a final concentration of 2% β -mercaptoethanol (284 mM). As an alternative, dithiothreitol (DTT or Cleland's reagent) may be used at a final concentration of 125 mM (19.3 mg/ml).

Note: For best results, do not store sample buffer with β -mercaptoethanol.

2. Dilute Sample

Dilute 1 part sample with 1 part sample buffer. More sample buffer can be added if necessary. A 1 part sample to 2 parts sample buffer dilution also works. Dry samples can be dissolved directly into the sample buffer.

Reference

Schagger H and von Jagow, Tricine-sodium dodecyl sulfate-polyacrylamide gel electrophoresis for the separation of proteins in the range from 1 to 100 kDa, Anal Biochem 166, 368-379 (1987)