

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

161-0734	10x Tris/Glycine , 1 L
161-0771	10x Tris/Glycine , 5 L

Native Sample Buffer

Catalog #

161-0738

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Dr., Hercules, CA 94547 USA

510-741-1000

4006027 Rev F

BIO-RAD

Introduction

Bio-Rad's native sample buffer is based on the method of Ornstein and Davis¹ (1964) with modifications specially formulated in our laboratory to improve band tightness. The use of native sample buffer ensures optimal band resolution when preparing proteins for polyacrylamide gel electrophoresis with Tris-glycine running buffer.

Specifications

Composition	62.5 mM Tris-HCl, pH 6.8 40% glycerol 0.01% Bromophenol Blue
Storage	4°C
Shelf life	1 year

Instructions for Use

Sample Dilution

Dilute 1 part sample with 1 part native 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

Ornstein L and Davis BJ, Disc Electrophoresis –1. Background and Theory, Ann NY Acad Sci 121, 321–349 (1964)