

Product Information

Catalog Number Product Description

Premixed Sample Buffers

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

Premixed Buffers

161-0733	10x TBE , 1 L
161-0770	10x TBE , 5 L
161-0732	10x Tris/Glycine/SDS , 1 L
161-0772	10x Tris/Glycine/SDS , 5 L
161-0734	10x Tris/Glycine , 1 L
161-0771	10x Tris/Glycine , 5 L
161-0744	10x Tris/Tricine/SDS , 1 L
161-0760	10x Tris/Tricine/SDS , 6 x 1 L

Bio-Rad Laboratories

2000 Alfred Nobel Dr., Hercules, CA 94547

510-741-1000

4106115 Rev B

TBE-Urea Sample Buffer

Catalog # 161-0768

The Bio-Rad logo consists of the words "BIO-RAD" in a bold, white, sans-serif font. The letters are contained within a black, rounded rectangular border that has a slight 3D effect, with a white highlight on the top edge.

Introduction

Bio-Rad's TBE-Urea Sample Buffer ensures optimal band resolution when preparing nucleic acids for denaturing polyacrylamide gel electrophoresis.

Specifications

<i>Composition</i>	<i>89mM Tris</i> <i>89mM Boric acid</i> <i>2mM EDTA, pH 8.0</i> <i>12% Ficoll</i> <i>0.01% Bromophenol Blue</i> <i>0.02% Xylene Cyanol FF</i> <i>7M Urea</i>
<i>Storage</i>	<i>4 °C</i>
<i>Shelf life</i>	<i>6 months</i>

Instructions For Use

Dilute 1 part sample with 1 part TBE-Urea Sample Buffer. Dry samples can be dissolved directly into the sample buffer. Denature samples by heating at 70 °C for 4 min prior to loading. Treatment at higher temperatures, up to 90 °C, may be required to completely denature some samples.