

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

Nucleic Acid Sample Loading Buffer, 5x

Catalog #
161-0767



Bio-Rad Laboratories

2000 Alfred Nobel Dr., Hercules, CA 94547

510-741-1000

4106116 Rev C

Introduction

Bio-Rad's Nucleic Acid Sample Loading Buffer ensures optimal band resolution when preparing nucleic acid samples for acrylamide and agarose gels. This sample buffer is formulated for use in both TBE and TAE buffer systems.

Specifications

<i>Composition</i>	<i>50 mM Tris-HCl, pH 8.0</i> <i>25% Glycerol</i> <i>5 mM EDTA</i> <i>0.2% Bromophenol Blue</i> <i>0.2% Xylene Cyanole FF</i>
<i>Storage</i>	<i>Ambient temperature</i>
<i>Shelf life</i>	<i>1 year</i>

Instructions For Use

Dilute 4-9 parts sample with 1 part Nucleic Acid Sample Buffer. Do not heat.

Reference

- 1 Maniatis, T., Fritsch, E.F., and Sambrook, J., *Molecular Cloning, A Laboratory Manual*, 160. Cold Spring Harbor Laboratory (1982)