

## Product Information

*Catalog Number Product Description*

### Premixed Sample Buffers

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

### Premixed Buffers

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