

## Product Information

*Catalog Number Product Description*

### Premixed Sample Buffers

161-0763	<b>IEF Sample Buffer</b> , 30 ml
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
161-0767	<b>Nucleic Acid Sample Buffer</b> , 5x, 10 ml
161-0768	<b>TBE-Urea Sample Buffer</b> , 30 ml

### Stain

161-0434	<b>IEF Staining Solution</b> , 1 L
----------	------------------------------------

### Premixed Buffers

161-0761	<b>10x Anode Buffer</b> , 250 ml
161-0762	<b>10x Cathode Buffer</b> , 250 ml

---

# Isoelectric Focusing (IEF) Gel Sample Buffer

## Catalog # 161-0763



### **Bio-Rad Laboratories**

2000 Alfred Nobel Dr., Hercules, CA 94547

510-741-1000

4006105 Rev C

## Introduction

Bio-Rad's IEF Gel Sample Buffer is composed of glycerol to allow sample loading of vertical IEF gels, such as Bio-Rad's Ready- and Criterion-brand gels.

## Specifications

*Sample Buffer  
Composition*                      *50% (w/w) Glycerol*

*Storage*                              *Ambient temperature*

*Shelf Life*                          *1 year*

## Instructions For Use

Dilute 9 parts sample with 1 part IEF Gel Sample Buffer. Dilutions resulting in a final glycerol concentration of less than 5% (i.e., a sample with more than 9 parts sample with 1 part sample buffer) are not recommended.