

# Product Information

*Catalog Number Product Description*

## **Premixed Sample Buffers**

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

## **Premixed Buffers**

161-0765	<b>10x Zymogram Renaturation Buffer, 125 ml</b>
161-0766	<b>10x Zymogram Development Buffer, 125 ml</b>
161-0732	<b>10x Tris/Glycine/SDS, 1 L</b>
161-0772	<b>10x Tris/Glycine/SDS, 5 L</b>

## **Bio-Rad Laboratories**

2000 Alfred Nobel Dr., Hercules, CA 94547

510-741-1000

4106111 Rev B

---

# Zymogram Sample Buffer

## Catalog #

# 161-0764



# Introduction

Bio-Rad's Zymogram Sample Buffer is based on the method of Laemmli.<sup>1</sup> The use of Zymogram Sample Buffer ensures optimal band resolution when preparing enzymes for Zymogram SDS-PAGE with Tris-Glycine-SDS running buffer.

# Specifications

<b>Composition</b>	<i>62.5 mM Tris-HCl, pH 6.8</i> <i>25% glycerol</i> <i>4% SDS</i> <i>0.01% Bromophenol Blue</i>
<b>Storage</b>	<i>Ambient temperature</i>
<b>Shelf life</b>	<i>1 year</i>

# Instructions For Use

Dilute 1 part sample with 2 parts Zymogram Sample Buffer. Dry samples can be dissolved directly into the sample buffer. Do not heat.

# Reference

<sup>1</sup> Laemmli, U.K., *Nature*, 227, 680 (1970).