

Product Information

Catalog # Description

Premixed Sample Buffers

- 161-0737 2x Laemmli Sample Buffer, 30 ml
- 161-0747 4x Laemmli Sample Buffer, 10 ml
- 161-0710 2-Mercaptoethanol, 25 ml
- 161-0738 Native Sample Buffer, 30 ml
- 161-0739 Tricine Sample Buffer, 30 ml
- 161-0767 5x Nucleic Acid Sample Buffer, 10 ml
- 161-0768 TBE-Urea Sample Buffer, 30 ml
- 161-0763 IEF Sample Buffer, 30 ml
- 161-0764 Zymogram Sample Buffer, 30 ml

Premixed Buffers

- 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

Bio-Rad Laboratories, Inc.

2000 Alfred Nobel Dr., Hercules, CA 94547 USA
510-741-1000

4006028 Rev J

2x Laemmli Sample Buffer

Catalog #161-0737

4x Laemmli Sample Buffer

Catalog #161-0747



BIO-RAD

Introduction

Bio-Rad's Laemmli sample buffers are based on the method of Laemmli (1970). The use of Laemmli sample buffers ensures optimal band resolution when preparing proteins for SDS-PAGE with Tris-glycine-SDS running buffer.

Formulations

2x Laemmli sample buffer	65.8 mM Tris-HCl, pH 6.8 26.3% (w/v) glycerol 2.1% SDS 0.01% bromophenol blue
4x Laemmli sample buffer	277.8 mM Tris-HCl, pH 6.8 44.4% (v/v) glycerol 4.4% LDS 0.02% bromophenol blue
Storage	Room temperature
Shelf life	2 years from date of manufacture

Instructions for Use

1. Add Reducing Agent

To obtain a final 1x concentration of 355 mM 2-mercaptoethanol

2x Laemmli sample buffer: Add 50 μ l of 2-mercaptoethanol per 950 μ l.

4x Laemmli sample buffer: Add 100 μ l of 2-mercaptoethanol per 900 μ l.

Alternatively, add dithiothreitol (DTT or Cleland's reagent) to a final 1x concentration of 50 mM.

Note: For best results, do not store sample buffer with 2-mercaptoethanol.

2. Dilute Sample

2x Laemmli sample buffer: Dilute 1 part sample with 1 part 2x Laemmli sample buffer.

4x Laemmli sample buffer: Dilute 3 parts sample with 1 part 4x Laemmli sample buffer. More sample buffer can be added if necessary.

Reference

Laemmli UK (1970). Cleavage of structural proteins during the assembly of the head of bacteriophage T4, Nature 227. 680–685.