

# electrophoresis standards

## POLYPEPTIDE STANDARDS

- SIZED FOR POLYPEPTIDE AND SMALL PROTEIN MOLECULAR WEIGHTS
- AVAILABLE PRESTAINED AND UNSTAINED
- STABLE FOR ONE YEAR

## SDS-PAGE and Kaleidoscope Standards

Bio-Rad's polypeptide standards are available as SDS-PAGE molecular weight standards or as prestained Kaleidoscope standards. These standards provide valuable information specifically for gel electrophoresis involving polypeptides and small molecular weight proteins.

### Determine Molecular Weights with SDS-PAGE Standards

The SDS-PAGE molecular weight standards are specifically designed for determining the molecular weights of peptides and small proteins on SDS-PAGE gels. Each tube contains six proteins, ranging from 1,423 to 26,625 daltons, which create a distinctive pattern of even intensity bands when separated on a Tris-Tricine gel. The standards produce excellent results when stained with a colloidal Coomassie<sup>®</sup> stain such as Coomassie G-250 dye or Bio-Safe<sup>™</sup> Coomassie stain.

### Monitor Runs with Kaleidoscope Standards

The Kaleidoscope Polypeptide Standards contain five individually colored proteins for easy identification of the protein bands. These standards are useful for assessing the transfer efficiency of a western blot, and for monitoring an electrophoretic gel during the run. The different colors make band identification very easy and serve as a quick reference to estimating the molecular weight range.

The molecular weight range of the five proteins is from 3,496 to 31,000 daltons. Each lot of Kaleidoscope Polypeptide Standards is calibrated using the SDS-PAGE polypeptide molecular weight standards and a calibration card is included with every lot.

### Both are Stable for One Year

SDS-PAGE and Kaleidoscope Polypeptide Standards are provided pre-blended in a stable storage buffer which allows the standards to be stored at -20 °C for up to 1 year without freeze/thaw degradation, eliminating the need to aliquot. Each vial is shipped with instructions and is guaranteed for 1 year when stored at -20 °C.

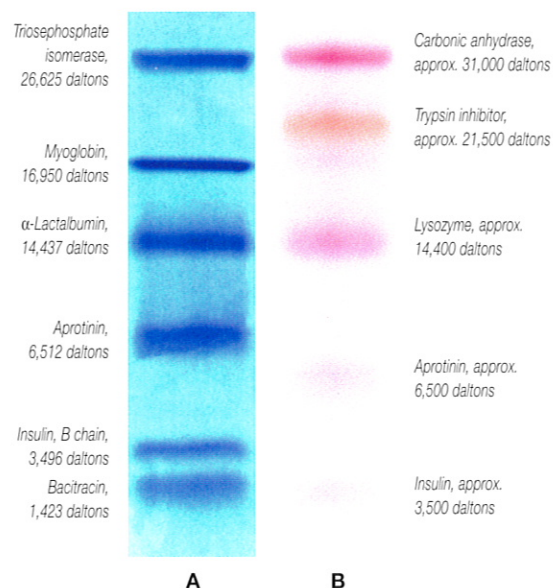


Fig. 1. (A) Polypeptide SDS-PAGE Standards run on a 16.5% Tris-Tricine Ready Gel and stained with Coomassie Blue G-250. (B) Kaleidoscope Polypeptide Standards.

**BIO-RAD**

## Specifications

	<b>SDS-PAGE</b>	<b>KALEIDOSCOPE</b>
<b>CONTENTS</b>	Rabbit triosephosphate isomerase Equine myoglobin Bovine $\alpha$ -lactalbumin Bovine aprotinin Bovine insulin, B chain, Oxidized bacitracin	Bovine carbonic anhydrase Soybean trypsin inhibitor Hen egg white lysozyme Bovine aprotinin Bovine insulin, B chain
<b>STORAGE BUFFER</b>	40% glycerol, 100 mM Tris-HCl pH 8.5, 4 mM EDTA, 3 mM NaN <sub>3</sub>	33% (v/v) glycerol, 0.5% SDS, 10 mM Tris, pH 7.0, 10 mM DTT, 2 mM EDTA, 0.01% NaN <sub>3</sub>
<b>VOLUME</b>	200 $\mu$ l	500 $\mu$ l
<b>STORAGE</b>	-20 °C	-20 °C
<b>SHIPPING CONDITIONS</b>	Room temperature	Room temperature
<b>SHELF LIFE</b>	1 year at -20 °C	1 year at -20 °C
<b>APPLICATIONS PER VIAL</b>	400	25–100
<b>RECOMMENDED GEL PERCENTAGES</b>	16.5% Tris-Tricine 10–20% Tris-Tricine	16.5% Tris-Tricine 10–20% Tris-Tricine

## Ordering Information

Catalog #	Description
161-0326	Polypeptide SDS-PAGE Standards, 200 $\mu$ l
161-0325	Kaleidoscope Polypeptide Standards, 500 $\mu$ l

*Coomassie is a trademark of ICI.*

**BIO-RAD**

**Bio-Rad  
Laboratories**

Life Science  
Group

Website [www.bio-rad.com](http://www.bio-rad.com) **U.S.** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Canada** (905) 712-2771  
**China** 86-10-62051850/51 **Denmark** 45 39 17 99 47 **Finland** 358 (0)9 804 2200 **France** 01 43 90 46 90 **Germany** 089 318 84-0  
**Hong Kong** 852-2789-3300 **India** (91-11) 461-0103 **Israel** 03 951 4127 **Italy** 39-02-216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460  
**Latin America** 305-894-5950 **Mexico** 514-2210 **The Netherlands** 0318-540666 **New Zealand** 64-9-4152280 **Norway** 22-74-18-70 **Russia** 7 095 979 98 00  
**Singapore** 65-2729877 **Spain** 34-91-661-7085 **Sweden** 46 (0)8-55 51 27 00 **Switzerland** 01-809 55 55 **United Kingdom** 0800-181134