

## Reference

1. Laemmli, U. K., *Nature*, **227**, 680 (1970).
2. Hames, B. D. and Rickwood, D., **Gel Electrophoresis of Proteins: A Practical Approach**, Second Edition, p. 17, Oxford University Press, New York (1990).

## Ordering Information

### Catalog

Number	Product Description
--------	---------------------

#### Molecular Weight Standards

161-0303	<b>SDS-PAGE Standards</b> , High, 200 $\mu$ l
161-0304	<b>SDS-PAGE Standards</b> , Low, 200 $\mu$ l
161-0317	<b>SDS-PAGE Standards</b> , Broad, 200 $\mu$ l
161-0314	<b>Silver Stain SDS-PAGE Standards</b> , Low, 200 $\mu$ l
161-0315	<b>Silver Stain SDS-PAGE Standards</b> , High, 200 $\mu$ l
161-0306	<b>Biotinylated SDS-PAGE Standards</b> , Low, 250 $\mu$ l
161-0311	<b>Biotinylated SDS-PAGE Standards</b> , High, 250 $\mu$ l
161-0319	<b>Biotinylated SDS-PAGE Standards</b> , Broad, 250 $\mu$ l
161-0320	<b>2-D SDS-PAGE Standards</b>
161-0326	<b>Polypeptide SDS-PAGE Standards</b> , 200 $\mu$ l

#### Prestained Standards

161-0305	<b>Prestained SDS-PAGE Standards</b> , Low, 500 $\mu$ l
161-0309	<b>Prestained SDS-PAGE Standards</b> , High, 500 $\mu$ l
161-0318	<b>Prestained SDS-PAGE Standards</b> , Broad, 500 $\mu$ l
161-0324	<b>Kaleidoscope Prestained Standards</b> , 500 $\mu$ l
161-0325	<b>Kaleidoscope Polypeptide Standards</b> , 500 $\mu$ l

#### IEF Standards

161-0310	<b>IEF Standards</b> , pI range 4.45-9.6, 250 $\mu$ l
----------	---

*Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules CA 94547*  
4006034 Rev C



# SDS-PAGE Molecular Weight Standards, High Range

Catalog Number  
**161-0303**

**Product shipped at room temperature.  
Store at -20 °C upon arrival.**

**BIO-RAD**

# Specifications

<b>Contents</b>	Approximately 400 µg of each protein blended to give bands of equal intensity on SDS polyacrylamide gels run according to Laemmli <sup>1</sup> and stained with Coomassie® blue R-250	
<b>Storage buffer</b>	50% glycerol, 300 mM NaCl, 10 mM Tris, 2 mM EDTA, 3 mM NaN <sub>3</sub>	
<b>Volume</b>	200 µl concentrated solution	
<b>Storage</b>	-20 °C	
<b>Shipping conditions</b>	Room temperature	
<b>Shelf life</b>	1 year at -20 °C	
<b>Applications per vial</b>	400 with Coomassie R-250	
<b>Recommended gel percentage*</b>	Low range	12.5%
	High range	7.5%
	Broad range	4-20 % gradient gels

**\*Note:** These standards can be run on other percentage gels, but all proteins may not be visible. Lower percentage gels may cause the low molecular weight proteins to migrate with or in front of the dye front. Higher percentage gels may prevent the high molecular weight proteins from separating.

## Protein Molecular Weights (daltons)

<b>Protein</b>	<b>Molecular Weight</b>	<b>High Range</b>	<b>Low Range</b>	<b>Broad Range</b>
Myosin	200,000	X		X
β-galactosidase	116,250	X		X
Phosphorylase b	97,400	X	X	X
Serum albumin	66,200	X	X	X
Ovalbumin	45,000	X	X	X
Carbonic anhydrase	31,000		X	X
Trypsin inhibitor	21,500		X	X
Lysozyme	14,400		X	X
Aprotinin	6,500			X

Coomassie is a trademark of ICI.

# Protocol

Dilute standards 1:20 in SDS Reducing Sample Buffer.\* Heat for 5 minutes at 95 °C. Cool and load 10 µl/well for full length gels (16-20 cm) or 5 µl/well for mini gels.

\* **SDS Reducing Sample Buffer** (prepare immediately before use)

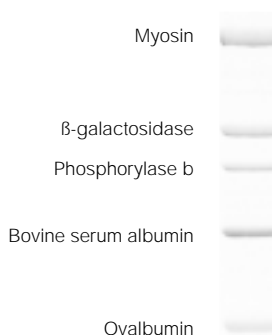
β-mercaptoethanol	25 µl
Stock Sample Buffer	475 µl
	500 µl

**Stock Sample Buffer** (store at room temperature)

Distilled water	4.8 ml
0.5 M Tris-HCl, pH 6.8	1.2 ml
Glycerol	1.0 ml
10% (w/v) SDS	2.0 ml
0.1% (w/v) Bromophenol blue	0.5 ml
	9.5 ml

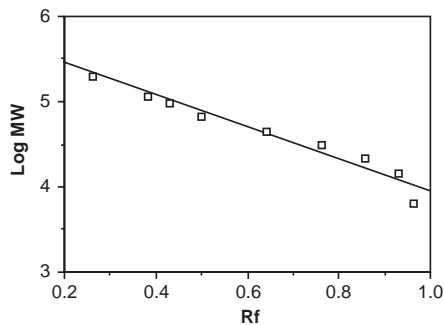
Use of Sample Buffer with insufficient or old β-mercaptoethanol may result in doublets at the soybean trypsin inhibitor and ovalbumin bands.

3



**Fig. 1. SDS polyacrylamide gels run in the Mini-PROTEAN® II cell according to the method of Laemmli.<sup>1</sup> High molecular weight standards run on a 7.5% SDS polyacrylamide gel, stained with Coomassie R-250.**

4



**Fig. 3. Curve generated by plotting the log of the molecular weight of the broad range standards vs. the relative mobility (Rf).**

$$R_f = \frac{\text{distance migrated by protein}}{\text{distance migrated by dye}}$$

The curve can be used to determine molecular weights of unknown proteins.<sup>2</sup>

5

## Protein References

Protein	Reference
<b>Rabbit skeletal muscle myosin</b>	Woods, E. F., Himmelfarb, S. and Harrington, W. F., <i>J. Biol. Chem.</i> , <b>238</b> , 2374 (1963).
<i>E. coli</i> β-galactosidase	Fowler, A. V. and Zabin, I., <i>Proc. Natl. Acad. Sci. USA</i> , <b>74</b> , 1507 (1977).
<b>Rabbit muscle phosphorylase b</b>	Titani, K., <i>et al.</i> , <i>Proc. Natl. Acad. Sci. USA</i> , Vol. <b>74</b> , 4762 (1977).
<b>Bovine serum albumin (BSA)</b>	Brown, J. R., <i>Fed. Proc.</i> , <b>34</b> , 591 (1975).
<b>Hen egg white ovalbumin</b>	Warner, R. C., "Egg Proteins," in: <b>The Proteins</b> , Vol. IIA, p. 435 (Neurath, H. and Bailey, K., eds.), Academic Press, New York (1954).
<b>Bovine carbonic anhydrase</b>	Davis, R. P., "Carbonic Anhydrase," in: <b>The Enzymes</b> , Vol V, p. 545. (Boyer, P. D., ed.) Academic Press, New York (1971).
<b>Soybean trypsin inhibitor</b>	Wu, Y. V. and Scheraga, H. A., <i>Biochemistry</i> , <b>1</b> , 698 (1962).
<b>Hen egg white lysozyme</b>	Jolles, P., <i>Angew. Chem Intl. Edit.</i> , <b>8</b> , 227 (1969).
<b>Bovine pancreatic trypsin inhibitor (Aprotinin)</b>	Kassell, B. and Laskowski, M., <i>Biochem. Biophys. Res. Comm.</i> , <b>20</b> , 463 (1965).

6