

Introduction

Bio-Rad's 2-D SDS-PAGE standards are formulated to provide a two-dimensional protein pattern with detection either by silver or Coomassie staining.

The standard proteins are defined by isoelectric point (pI) and molecular weight. Added to the sample, they are used to determine pIs and molecular weights of sample proteins or to serve as a marker for 2-D gel matching. The characteristic pattern of main and minor spots of each standard protein makes them easy to identify among sample spots.

In addition, the standards can be used alone as an internal control for reproducibility of 2-D experiments or to test ampholyte mixtures.

The most striking advantage of the standards is that they allow valid comparison of 2-D electrophoretic patterns between different laboratories.

2-D SDS-PAGE Standards, Specifications

pI range	4.5 to 8.5
Molecular weight range	17,500 to 76,000
Contents	Hen egg white conalbumin type 1 Bovine serum albumin (BSA) Bovine muscle actin Rabbit muscle glyceraldehyde 3-phosphate dehydrogenase (GAPDH) Bovine carbonic anhydrase Soybean trypsin inhibitor Equine myoglobin 9 M urea 5% 2-mercaptoethanol 2% Bio-Lyte® 5/7 ampholyte Distilled water
Volume	500 µl solution, ready to use
Shipping	4°C
Storage	-20°C Aliquoting is recommended prior to freezing. Frequent freeze/thaw procedures should be avoided. Aliquots can be stored at 4°C for about one week
Shelf life	One year
Applications per vial	Important! Always vortex the 2-D SDS-PAGE standards after thawing or refrigerating to resuspend the urea 200 (2.5 µl per mini gel) using Coomassie stain 1,000 (0.5–2.5 µl per mini gel) using silver stain Double these amounts for full-length gels (16–20 cm)

Ordering Information

Catalog #	Description
161-0320	2-D SDS-PAGE Standards
161-0303	SDS-PAGE Standards, high range
161-0304	SDS-PAGE Standards, low range
161-0363	Precision Plus Protein™ Unstained Standards
161-0373	Precision Plus Protein All Blue Standards
161-0374	Precision Plus Protein Dual Color Standards

Instruction Manual

2-D SDS-PAGE Standards

BIO RAD

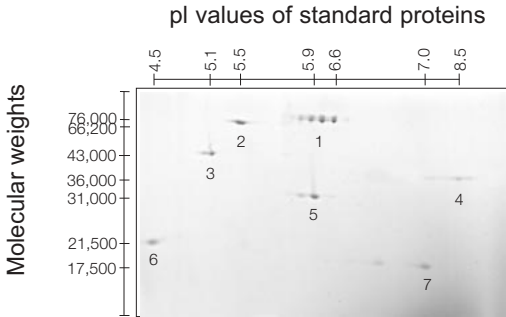
Bio-Rad Laboratories, Inc.

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

4110064 Rev A

BIO RAD

Two-Dimensional Electrophoretic Protein Pattern



Protein spots

- 1 Conalbumin
- 2 BSA
- 3 Actin
- 4 GAPDH
- 5 Carbonic anhydrase
- 6 Trypsin inhibitor
- 7 Myoglobin

Two-dimensional electrophoretic protein pattern of 2-D SDS-PAGE standards separated in the Mini-PROTEAN® II cell. 2.5 µl of the standards were run according to the method of Klose (Klose J, Humangenetik 26, 231–243, 1975; Klose K and Feller M, Electrophoresis 2, 12–24, 1981; Jungblut PR and Seifert R, J Biochem Biophys Methods 21, 47–58, 1990)

Protein Molecular Weights and Isoelectric Points (pI)

Protein	Molecular Weight	pI	Reference
Hen egg white conalbumin type 1	76,000	6.0	Cavorta P, et al., Experientia 34, 849–850 (1978)
		6.3	
		6.6	
Bovine serum albumin (BSA)	66,200	5.4	Brown JR, Fed Proc 34, 591 (1975)
		5.6	
Bovine muscle actin	43,000	5.0	The running pI value of muscle actin has been empirically determined.
		5.1	
Rabbit muscle GAPDH	36,000	8.3	Smith CM and Velick SF, J Biol Chem 247, 273–284, (1972)
		8.5	

Protein	Molecular Weight	pI	Reference
Bovine carbonic anhydrase	31,000	5.9	Ashworth RB, et al., Arch Biochem Biophys 142, 122–131 (1971) Jonsson M and Petterson E, Acta Chem Scand, 22, 712–713 (1966) Davis RP, Carbonic Anhydrase, p. 545–564 in The Enzymes, Vol V, 2nd edn (Boyer PD, ed), Academic Press, New York (1971)
		6.0	
Soybean trypsin Inhibitor	21,500	4.5	Catsimpoulas N and Leuthner E, Anal Biochem 31, 437–447 (1969) Wu YV and Scheraga HG, Biochemistry 1, 698–705 (1962)
Equine myoglobin	17,500	7.0	Salaman MR and Williamson AR, Biochem J 122, 93–99 (1971)