

Ordering Information

Catalog

Number Product Description

Specialty Standards

161-0310 **IEF Standards**, pl range 4.45-9.6, 250 μ l

161-0320 **2-D SDS-PAGE Standards**, 500 μ l

Prestained Standards

161-0324 **Kaleidoscope Prestained Standards**, 500 μ l

161-0325 **Kaleidoscope Polypeptide Standards**, 500 μ l

161-0305 **Prestained SDS-PAGE Standards**,
low range, 500 μ l

161-0309 **Prestained SDS-PAGE Standards**,
high range, 500 μ l

161-0318 **Prestained SDS-PAGE Standards**,
broad range, 500 μ l

SDS-PAGE Molecular Weight Standards

161-0304 **SDS-PAGE Standards**, low range, 200 μ l

161-0303 **SDS-PAGE Standards**, high range, 200 μ l

161-0317 **SDS-PAGE Standards**, broad range, 200 μ l

161-0326 **Polypeptide Standards**, 200 μ l

161-0314 **Silver Stain Standards**, low range, 200 μ l

161-0315 **Silver Stain Standards**, high range, 200 μ l

161-0306 **Biotinylated Standards**, low range, 250 μ l

161-0311 **Biotinylated Standards**, high range, 250 μ l

161-0319 **Biotinylated Standards**, broad range, 250 μ l



IEF Standards Instruction Manual

Catalog Number 161-0310
Broad Range pl 4.45-9.6

Product shipped at room temperature.

For research use only.

Store at -20 °C.

BIO-RAD

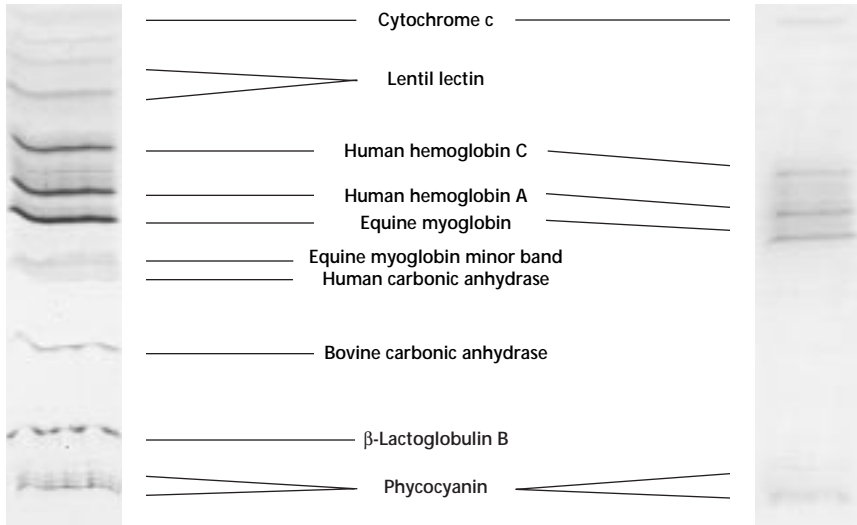


Fig. 1. IEF Standards for accurate pI calibration of native IEF gels. A. 5 μ l of the standards were run on a horizontal polyacrylamide IEF gel and stained with Coomassie blue R-250 dye and crocein scarlet. B. 5 μ l of the standards were run on a horizontal polyacrylamide IEF gel and left unstained.

Introduction

Bio-Rad's IEF Standards, a mixture of nine natural proteins with isoelectric points ranging from 4.45 to 9.6, permit dependable and reproducible *pI* calibration in IEF gels. The mixture is provided in a stable aqueous solution. No reconstitution or dilution is required prior to use. IEF Standards are intended for use in analytical acrylamide or agarose isoelectric focusing gels. Five of the nine proteins are naturally colored to provide continuous monitoring of the focusing process.

3

Constituent Proteins

	Color	<i>pI</i>	MW
Phycocyanin (3 bands)	Blue	4.45, 4.65, 4.75	232,000
β-Lactoglobulin B		5.1	18,400
Bovine carbonic anhydrase		6.0	31,000
Human carbonic anhydrase		6.5	28,000
Equine myoglobin (2 bands)	Brown	6.8, 7.0	17,500
Human hemoglobin A	Red	7.1	64,500
Human hemoglobin C	Red	7.5	64,500
Lentil lectin (3 bands)		7.80, 8.00, 8.20	49,000
Cytochrome c	Red	9.6	12,200

Note: The *pI* values given here were determined by direct measurement with a surface pH electrode.

5

Specifications

Contents	Approximately 16.5 mg total protein in 50% glycerol with 0.02% sodium azide
Volume	250 µl
Storage	-20 °C*
Shelf life	1 year at -20 °C
Recommended sample volume	5 µl for Coomassie blue staining (0.5 µl for silver stain)
Applications per vial	50 (500 for silver stain)

* IEF Standards will be shipped at room temperature. They are stable for transportation at this temperature. Store at -20 °C upon arrival. Do not store at -70 °C.

Note: IEF Standards are not recommended for 2-D electrophoresis applications. The proteins are not naturally colored when denatured by urea and subsequent electrophoresis will produce different results than those shown in Figure 1.

4

Instructions for Use

IEF Standards can be applied directly to IEF gels without any prior treatment. Use 5 µl for gels which are to be stained with Coomassie blue, and for silver staining, use 0.5 µl or 5 µl of a 1:10 dilution of IEF Standards, using deionized, distilled water as a diluent.

6