

---

# ProteinChip® Protein Calibrant Kit

Catalog #C10-00001

## Kit Contents

Protein Standard	Amount/vial	# vials	Molecular Weight
Insulin, bovine	1 nmol	2	5,733.6
Ubiquitin, bovine	1 nmol	2	8,564.8
Cytochrome C, bovine	1 nmol	2	12,230.9
Superoxide dismutase, bovine	1 nmol	2	15,591.4
Myoglobin, equine	1 nmol	2	16,951.5
β-lactoglobulin A, bovine	1 nmol	2	18,363.3
Horseradish peroxidase	2 nmol	2	43,240.0
Serum albumin, bovine	2 nmol	2	66,410.0
Conalbumin, chicken	2 nmol	2	77,490.0
IgG, bovine	5 nmol	2	147,300.0

## Preparation (allow 30 min)

1. Briefly centrifuge vials containing dried protein standards to ensure proteins are at the bottom of the vials.
2. Add diluent to protein standard vials.
  - For insulin protein standard, aliquot 100 µl 10 mM HCl into vial
  - For all other standards, aliquot 100 µl deionized water into each vial
3. Cap and vortex each vial for several seconds; let stand 30 min, then vortex again to dissolve protein.

## Storage

Store all vials containing protein, whether dry or dissolved, at -20°C.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 4BIORAD **Australia** 61 02 9914 2800  
**Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400 **Canada** 905 712 2771  
**China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00  
**Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 318 84 0  
**Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800  
**India** 91 124 4029300 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 5811 6270  
**Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 0508 805 500 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99  
**Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188  
**South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700  
**Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000