
ProteinChip® Peptide Calibrant Kit

Catalog #C10-00002

Kit Contents

Peptide Standard	Amount/vial	# vials	Molecular Weight
Angiotensin I, human	1 nmol	2	1,296.5
[Glu ¹]-fibrinopeptide B	1 nmol	2	1,570.6
Dynorphin A (209–225), porcine	1 nmol	2	2,147.5
ACTH (1–24), human	1 nmol	2	2,933.5
β-endorphin (61–91), human	1 nmol	2	3,465.0
Insulin, bovine	1 nmol	2	5,733.6
Ubiquitin, bovine	1 nmol	2	8,564.8

Preparation (allow 30 min)

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

Storage

Store all vials containing peptides, whether dry or dissolved, at –20°C.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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