

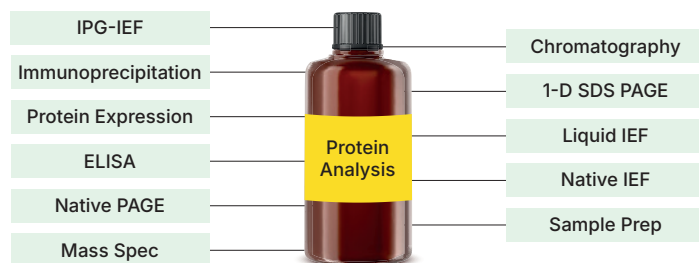
Reagent Compatibility Chart for Bio-Rad Protein Assays

Reagent Compatibility

The following reagents were tested for their compatibility with Bio-Rad protein assays: Quick Start Bradford Protein Assay, Bio-Rad Protein Assay, DC Protein Assay, and RC DC Protein Assay. The concentrations represent maximum concentrations for standard assay.

	Quick Start Bradford Protein Assay	Bio-Rad Protein Assay	DC Protein Assay	RC DC Protein Assay
Acetate	—	0.6 M	—	—
Acetone	10%	✓	—	—
Acetonitrile	10%	—	—	—
Acidic pH	—	✓	—	—
Adenosine	—	0.001 M	—	—
Amino acids	—	—	✓	—
Ammonium sulfate	1 M	1 M	0.5 M	—
Ampholytes pH 3–10	0.5%	0.5%	—	—
ASB-14	0.03%	—	—	—
Ascorbic acid	0.05 M	—	—	—
ATP	—	0.001 M	—	—
Barbital	—	✓	—	—
BES	—	2.5 M	—	—
Bis-Tris, pH 6.5	0.2 M	—	—	—
Boric acid	—	✓	—	—
Brij-35	—	—	1%	—
C ₁₂ E ₈	—	—	0.2%	—
Cacodylate-Tris	—	0.1 M	—	—
Calcium chloride	0.04 M	—	0.05 M	—
CDTA	—	0.05 M	—	—
CHAPS	10%	—	1%	2%
CHAPSO	10%	—	1%	1%
Citrate	—	0.05 M	—	—
Deoxycholate	—	0.1%	—	—
Deoxycholic acid	0.2%	—	—	—
Dithioerythritol (DTE)	0.01 M	—	—	—
Dithiothreitol (DTT)	0.01 M	1 M	0.001 M	0.1 M
DMSO	5%	—	—	—

Where to Use a Protein Assay



	Quick Start Bradford Protein Assay	Bio-Rad Protein Assay	DC Protein Assay	RC DC Protein Assay
DNA	—	1 mg/ml	—	—
Eagle's MEM	✓	✓	—	—
Earle's salt solution	✓	✓	—	—
EDTA	0.2 M	0.1 M	0.025 M	0.1 M
EGTA	0.2 M	0.05 M	—	—
Ethanol	10%	✓	—	—
Formic acid	—	1.0 M	—	—
Fructose	—	✓	—	—
Glucose	20%	✓	—	—
Glutathione	—	✓	—	—
Glycerol	5%	99%	—	—
Glycine	0.1 M	0.1 M	—	—
Guanidine HCl	2 M	✓	0.4 M	—
Hank's salt solution	✓	✓	—	—
HCl	0.1 M	—	0.5 M	—
HEPES	0.1 M	0.1 M	—	—
Imidazole	0.2 M	—	—	0.5 M
Laemmli buffer	—	—	—	✓
Magnesium chloride	1 M	1 M	—	—
Malic acid	—	0.2 M	—	—
2-Mercaptoethanol	1 M	1 M	×	5%
MES	0.1 M	0.7 M	—	—
Methanol	10%	✓	—	—
Modified Dulbecco's PBS	✓	—	—	—
MOPS	0.1 M	0.2 M	—	—
NAD	0.002 M	0.001 M	—	—
NaSCN	—	3 M	—	—
Nonidet P-40	0.25%	—	2%	—

✓, compatible; —, not tested; ×, not compatible; DC, detergent compatible; RC DC, reducing agent and detergent compatible.



	Quick Start Bradford Protein Assay	Bio-Rad Protein Assay	DC Protein Assay	RC DC Protein Assay
Octyl β-glucoside	0.5%	—	1%	—
Octyl β-thioglucopyranoside	1%	—	—	—
PBS	✓	—	—	—
Peptone	—	✓	—	—
Phenol Red	0.5 mg/ml	—	—	—
Phenol	—	5%	—	—
Phosphate	—	1 M	—	—
PIPES	0.2 M	0.5 M	—	—
PMSF	0.002 M	—	—	—
Polyadenylic acid	—	0.001 M	—	—
Polypeptides (MW <3000)	—	✓	—	—
Potassium chloride	2 M	1 M	—	—
Potassium phosphate	0.5 M	—	—	—
Pyrophosphate	—	0.2 M	—	—
rRNA	—	0.25 mg/ml	—	—
SB 3-10	0.1%	—	—	—
SDS	0.03%	0.1%	10%	—
Sodium acetate, pH 4.8	0.2 M	—	—	—
Sodium azide	0.5%	—	0.05%	—
Sodium bicarbonate	0.2 M	—	—	—
Sodium carbonate	0.1 M	—	—	—
Sodium chloride	2.5 M	5 M	—	—
Sodium citrate, pH 4.8 or 6.4	0.2 M	—	—	—

	Quick Start Bradford Protein Assay	Bio-Rad Protein Assay	DC Protein Assay	RC DC Protein Assay
Sodium hydroxide	0.1 M	—	0.5 M	2.5 M
Sodium phosphate	0.5 M	✓	—	—
Streptomycin sulfate	—	20%	—	—
Sucrose	10%	—	—	—
TBP	0.005 M	—	—	0.002 M
TBS	0.5x	—	—	—
TCEP	0.02 M	—	—	—
Thesit	—	—	1%	—
Thiourea	1 M	—	—	—
Thymidine	—	0.001 M	—	—
Total RNA	—	0.30 mg/ml	—	—
Tricine	0.05 M	✓	—	—
Triethanolamine, pH 7.8	0.05 M	—	—	—
Tris, pH 8	1 M	2 M	0.1 M	0.5 M
Tris/glycine	1x	—	—	—
Tris/glycine/SDS	0.5x	—	—	—
Triton X-100	0.05%	0.1%	1%	2%
tRNA	—	0.4 mg/ml	—	—
Tween 20	0.01%	—	1%	2%
Tyrosine	—	0.001 M	—	—
Urea	4 M	6 M	4 M	—
Vitamins	—	✓	—	—

✓, compatible; —, not tested; ×, not compatible; DC, detergent compatible; RC DC, reducing agent and detergent compatible.

Ordering Information

Catalog #	Description
5000001	Bio-Rad Protein Assay Kit I, includes dye concentrate and bovine γ-globulin standard
5000002	Bio-Rad Protein Assay Kit II, includes dye concentrate and bovine serum albumin standard
5000201	Quick Start Bradford Protein Assay Kit 1, includes 1 L 1x dye reagent, BSA standard (5 x 2 mg/ml)
5000202	Quick Start Bradford Protein Assay Kit 2, includes 1 L 1x dye reagent, BSA standard set (2 sets of 7 standards, 2 ml of 0.125–2.0 mg/ml)
5000203	Quick Start Bradford Protein Assay Kit 3, includes 1 L 1x dye reagent, bovine γ-globulin standard (5 x 2 mg/ml)
5000204	Quick Start Bradford Protein Assay Kit 4, includes 1 L 1x dye reagent, bovine γ-globulin standard set (2 sets of 7 standards, 2 ml of 0.125–2.0 mg/ml)

Catalog #	Description
5000111	DC Protein Assay Kit I, includes DC protein assay reagents package and bovine γ-globulin standard
5000112	DC Protein Assay Kit II, includes DC protein assay reagents package and bovine serum albumin standard
5000121	RC DC Protein Assay Kit I, includes RC reagents package, DC protein assay reagents package, bovine γ-globulin standard
5000122	RC DC Protein Assay Kit II, includes RC reagents package, DC protein assay reagents package, bovine serum albumin standard

Visit [bio-rad.com/proteinassaykits](https://www.bio-rad.com/proteinassaykits) for more information.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc.
All trademarks used herein are the property of their respective owner.
© 2025 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science Group

Website [bio-rad.com](https://www.bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399 Canada 1 800 361 1808 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23 France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050 Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670 The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23 Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23 Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23