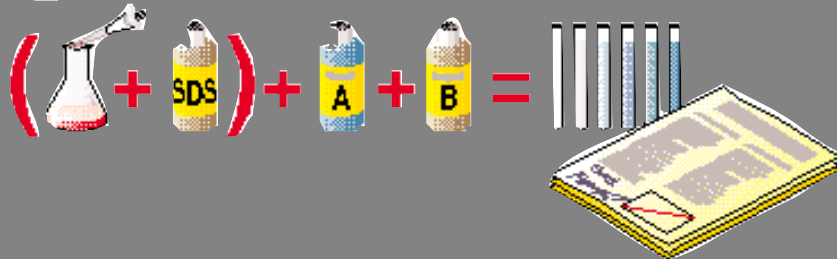
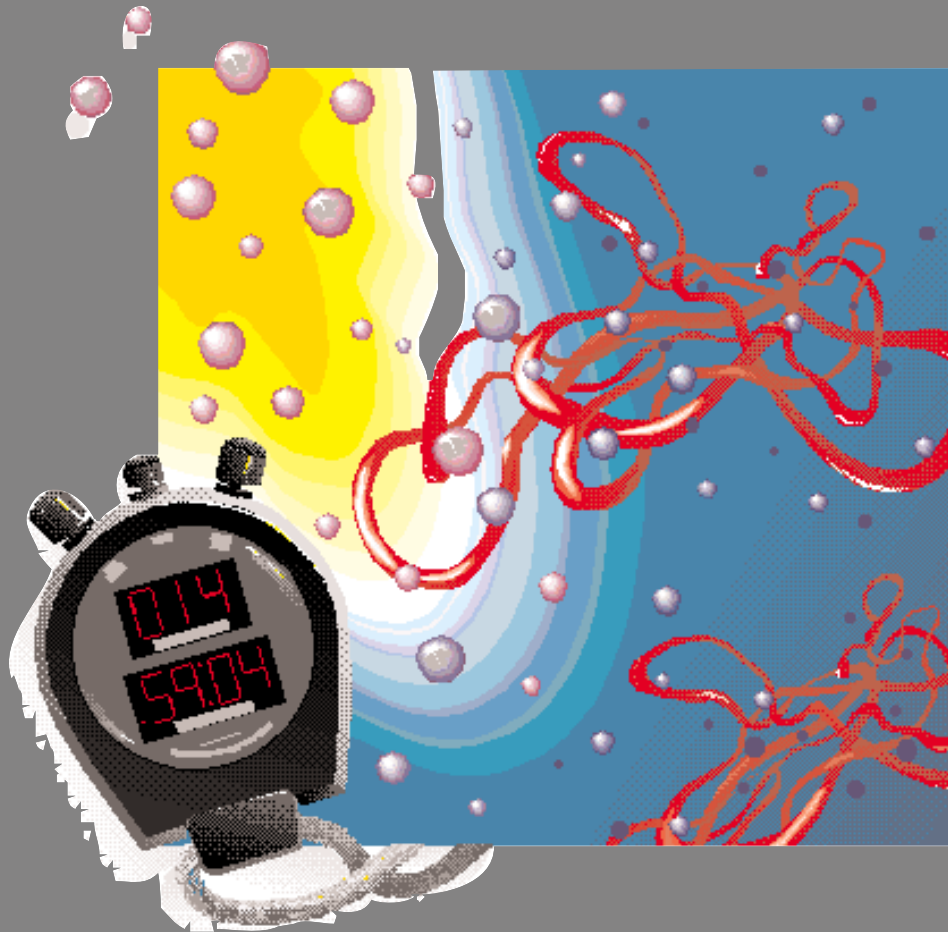




DC Protein Assay



BIO-RAD



The Detergent- Compatible (DC) Assay for Protein

Bio-Rad's DC Protein Assay is specially suited for samples solubilized in detergent. A proprietary modification of the traditional Lowry method, the DC Protein Assay measures detergent-solubilized proteins quickly and reproducibly following a single 15 minute incubation.

Increased color stability

The color will not change more than 5% in 1 hour, or 10% in 2 hours. This means that, in spite of inevitable interruptions, your standard curve is good from the first to the last sample.

Simplified, detergent-compatible protein assay

After 15 minutes at room temperature, samples are ready to be evaluated. No timed, multi-step reagent additions or incubations in the water bath are required.

The chemistry is compatible with up to 1% ionic or nonionic detergent in addition to other common reagents such as EDTA and Tris (see Reagent Compatibility chart).

Convenience in a kit

The Bio-Rad DC Protein Assay reagents provide all the necessary reagents for 500 standard assays or 10,000 microplate assays. The reagents may be purchased with an IgG standard, a BSA standard, or by themselves in the DC Protein Assay Reagents Package.

Free technical bulletin available

Simply call your local Bio-Rad representative (1-800-4BIORAD in the U.S.) and

request bulletin 1069. It describes both the Bio-Rad (Bradford) Protein Assay and the DC Protein Assay.

Reagent Compatibility

10% SDS	1% CHAPS*
2% NP-40	1% Triton® X-100
1% CHAPSO	1% octyl glucoside
1% Thesit®	1% Tween® 20
1% Brij®-35	0.2% C ₁₂ E ₈ *
0.1 M Tris, pH 8.0	0.5 M NaOH
0.5 M HCl	0.5 M (NH ₄) ₂ SO ₄
0.025 M EDTA	0.05 M CaCl ₂
0.4 M Guanidine HCl	4 M Urea
0.05% Sodium azide	1% Triton X-114
1% deoxycholate	1 mM DTT
0.3% digitonin	

The listed compounds were tested and found to be compatible with the Bio-Rad DC Protein Assay. In some cases, the presence of one or more of these substances will change the response of the protein to the assay reagents; therefore, the standard should always be prepared in the same buffer as the sample.

	Bio-Rad Protein Assay	Bio-Rad DC Protein Assay
Standard assay sample size	100 µl	100 µl
Microplate assay sample size	10 µl	5 µl
Microassay sample size	800 µl	200 µl
Linear range for standard assay	0.2–1.4 mg/ml	0.1–2.0 mg/ml
Linear range for microassay	1.25–25 µg/ml	5–250 µg/ml
Minimum incubation time	5 minutes	15 minutes
Unique compatibility	Reducing agents	Detergents, basic protein buffers
Standard assays per kit	450	500
Microplate assays per kit	2,250	10,000
Adapted from method by	Bradford	Lowry

Ordering Information

Catalog Number Description

For samples solubilized in detergent:

500-0111	DC Protein Assay Kit I , includes reagents and IgG standard
500-0112	DC Protein Assay Kit II , includes reagents and BSA standard
500-0116	DC Protein Assay Reagents Package , includes reagents without a standard

For determinations by method of Bradford:

500-0001	Protein Assay Kit I , includes dye concentrate and IgG standard
500-0002	Protein Assay Kit II , includes dye concentrate and BSA standard
500-0006	Dye Reagent Concentrate , includes reagent without a standard

†Brij and Tween are registered trademarks of Atlas Chemical. Thesit is a registered trademark of Desitin Arzneimittel GmbH. Triton is a registered trademark of Rohm and Haas.

*octaethyleneglycol dodecyl ether

BIO-RAD

Bio-Rad
Laboratories

Life Science
Group

U.S. (800) 4BIORAD • California Ph. (510) 741-1000 • New York Ph. (516) 756-2575 • Australia Ph. 02-805-5000 • Austria Ph. (1)-877 89 01 • Belgium Ph. 09-385 55 11 • Canada Ph. (905) 712-2771 • China Ph. (01) 2051873 • France Ph. (1) 49 60 68 34 • Germany Ph. 089 31884-0 • Italy Ph. 02-21609 1 • Japan Ph. 03-3534-7665 • Hong Kong Ph. 7893300 • The Netherlands Ph. 08385-40666 • New Zealand Ph. 09-443 3099 • Singapore Ph. (65) 4432529 • Spain Ph. (91) 661 70 85 • Sweden Ph. 46 (0) 8-735 83 00 • Switzerland Ph. 01-810 16 77 • United Kingdom Ph. 0800 181134