



Bio-Safe™ Coomassie Stain

Eliminate the Use of
Hazardous Chemicals.

- No solvent waste disposal costs
- No hazardous shipping charges
- No special handling
- No fume hood requirement

For more information, contact your local Bio-Rad representative
or visit us on the Web at discover.bio-rad.com



BIO-RAD

Bio-Safe Coomassie Stain



Bio-Safe colloidal Coomassie Blue G-250 stain from Bio-Rad is a premixed solution that is nonhazardous and does not require the use of methanol and acetic acid for destaining. Bio-Safe stain produces blue bands on a clear background, with sensitivity to 8 ng of protein.

Fast, Simple, Sensitive, and Convenient

- Fast (60–120 min from start to finish)
- Premixed ready-to-use solution in both 1 L and 5 L sizes (20–50 ml per mini gel)
- Detection range of 8–29 ng per band
- Bands/spots appear and can be viewed while gel is in the stain
- No gel shrinkage due to methanol
- Water wash prevents over-destaining
- Compatible with an alcohol fixative step, if desired
- Mass spectrometry compatible

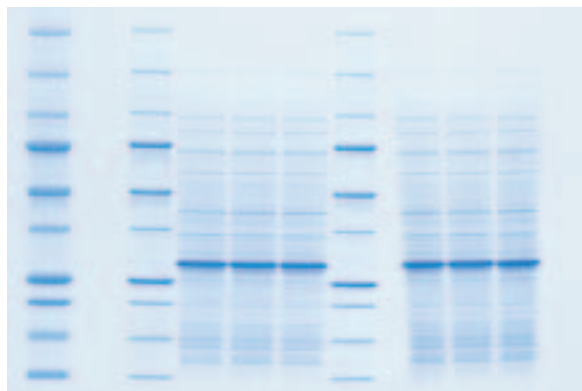
Broad Applications

- SDS-PAGE
- Native PAGE
- Peptide analysis (Tricine gels with fixation step)
- 2-D electrophoresis

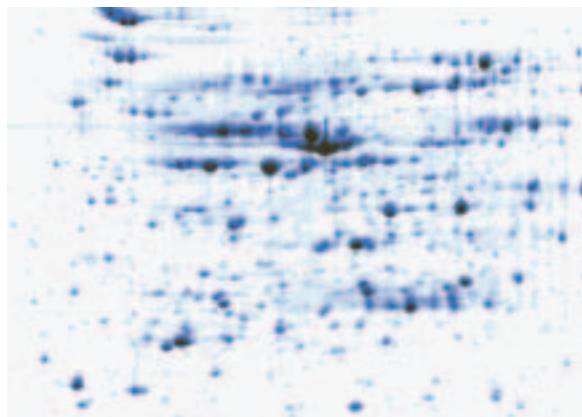
Ordering Information

Catalog #	Description
161-0786	Bio-Safe Coomassie Stain, 1 L bottle, stains 20 mini gels
161-0787	Bio-Safe Coomassie Stain, 5 L cube, stains 100 mini gels

Coomassie is a trademark of BASF Aktiengesellschaft.



1-D gel stained with Bio-Safe Coomassie stain. A Criterion™ 4–20% gel was washed in water for 20 min, stained with Bio-Safe Coomassie stain for 60 min, then destained in water for 60 min.



2-D gel stained with Bio-Safe Coomassie stain. First dimension: 300 µg of *E. coli* lysate treated with RNase and DNase in rehydration buffer containing 0.1% Bio-Lyte® ampholytes was loaded onto a 17 cm ReadyStrip™ IPG strip (pH 4–7) and focused 92,000 V-hr on the PROTEAN® IEF cell. Second dimension: The strip was loaded onto a 20 x 20.5 cm PROTEAN® Plus 8–16% gel, and run at 200 V constant for 6 hr.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Brazil** 55 21 2527 3454 **Canada** (905) 712-2771 **China** (86-21) 63052255 **Czech Republic** + 420 2 41 43 05 32 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 318 84-0 **Hong Kong** 852-2789-3300 **Hungary** 36 1 455 8800 **India** (91-124)-6398112/113/114, 6450092/93 **Israel** 03 951 4127 **Italy** 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 55-52-00-05-20 **The Netherlands** 0318-540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** + 48 22 331 99 99 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-6415 3188 **South Africa** 00 27 11 4428508 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717-9555 **Taiwan** (8862) 2578-7189/2578-7241 **United Kingdom** 020 8328 2000