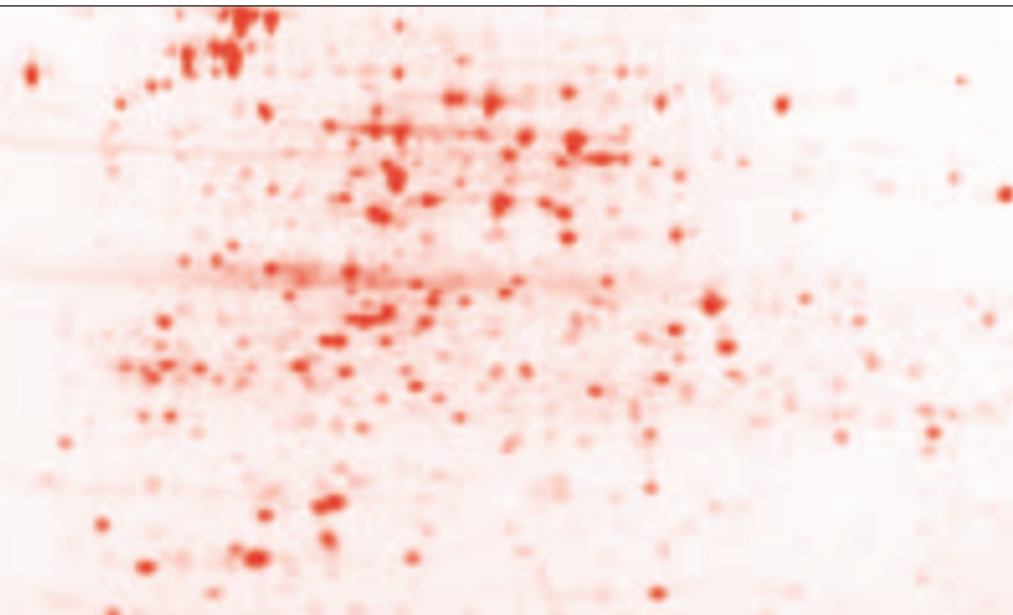


SYPRO Ruby Protein Gel Stain



linearity



sensitivity



versatility



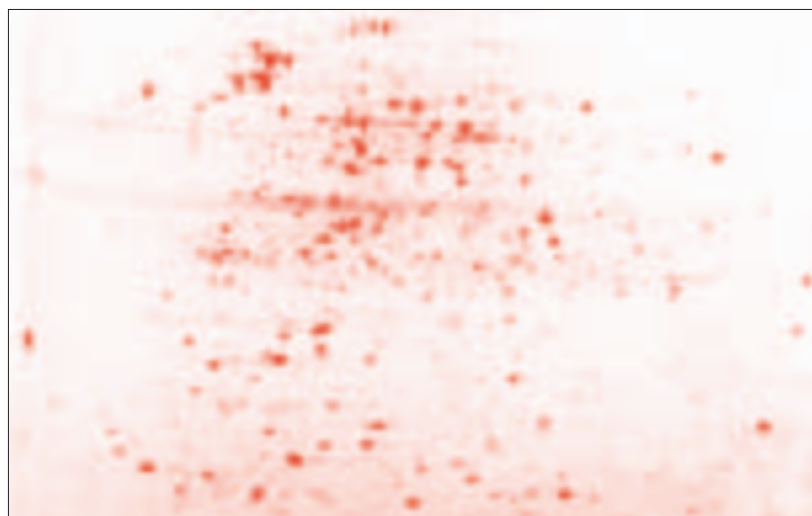
simplicity



The Answer to Your **2-D Staining** Needs

SYPRO Ruby protein gel stain is a fluorescent stain for polyacrylamide gels that extends the range of staining over 3 orders of magnitude, is adaptable to high-throughput applications, and is compatible with downstream microanalysis (mass spectrometry and Edman-based sequencing). The protocol provides walk-away convenience and nanogram-level sensitivity with no overstaining.

BIO-RAD



E. coli extract (100 µg) was focused on a ReadyStrip™ IPG strip (pH 3–10), separated on a 12% Criterion™ Tris-HCl gel, then stained with SYPRO Ruby protein gel stain.

SYPRO Ruby Protein Gel Stain

Linearity

- Reproducible protein quantitation over 3 orders of magnitude

Simplicity

- Single-reagent, ready-to-use formula
- Flexible staining time (3 hr to overnight) with no risk of overstaining

Versatility

- Compatible with subsequent analyses (mass spectrometry, blotting, and Edman-based sequencing)
- Excitable by blue or UV wavelengths
- Provides maximum sensitivity and image resolution with laser scanners

Sensitivity

- Sensitivity equivalent to silver stain (1–10 ng)
- Specific for proteins; no staining of nucleic acids

Bibliography

Berggren K et al., Background-free, high sensitivity staining of proteins in one- and two-dimensional sodium dodecyl sulfate-polyacrylamide gels using a luminescent ruthenium complex, *Electrophoresis* 21, 2509–2521 (2000)

Ordering Information

Catalog #	Description
170-3126	SYPRO Ruby Protein Gel Stain, 200 ml
170-3125	SYPRO Ruby Protein Gel Stain, 1 L
170-3138	SYPRO Ruby Protein Gel Stain, 5 L cube
170-3127	SYPRO Ruby Protein Blot Stain, 200 ml

SYPRO is a trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell SYPRO products for research use only, under US patent 5,616,502.

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