

## Order Today

Enjoy the benefits of using SYPRO Orange. Order today, or contact your local Bio-Rad office to find out more about this unique stain.

### Catalog No. Product Description

170-3120	<b>SYPRO Orange Concentrate,</b> 500 µl, sufficient for 50–100 minigels
161-0332	<b>SDS-PAGE Standards for SYPRO Orange Stain,</b> broad range, 6.5–200 kd, 200 µl (other size ranges available)

This product is offered for research purposes only. Coomassie is a registered trademark of ICI Organics, Inc. Polaroid is a registered trademark of Polaroid Corp. SYPRO is a registered trademark of, and the dye is manufactured by, Molecular Probes, Inc. Patents pending.



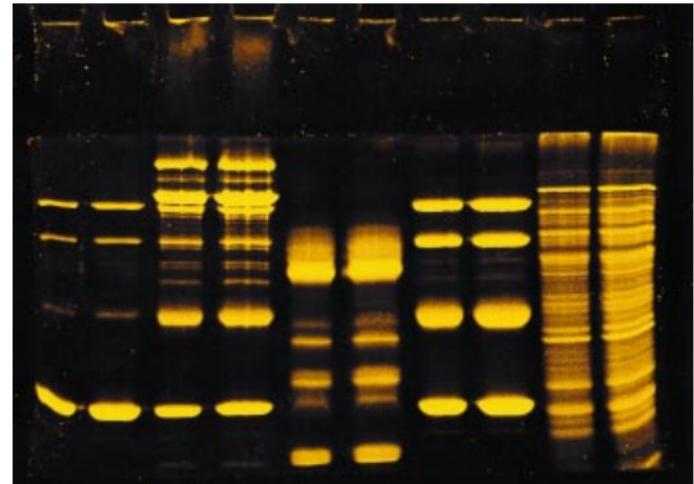
**Bio-Rad**  
Laboratories

**Molecular  
Bioscience Group**

U.S. (800) 4BIORAD California (510) 741-1000 Australia 02-9914-2800 Austria (1) 877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771 China (01) 2046622 Denmark 39 17 9947 Finland 90 804 2200 France (1) 49 60 68 34 Germany 089 31884-0 India 91-11-461-0103 Italy 02-21609 1 Japan 03-5811-6270 Hong Kong 7893300 The Netherlands 0318-540666 New Zealand 09-443 3099 Singapore (65) 272-9877 Spain (91) 661 70 85 Sweden 46 (0) 735 83 00 Switzerland 01-809 55 55 United Kingdom 0800 181134

# New! SYPRO® Orange

## High-Sensitivity Fluorescent Protein Gel Stain



One 30-minute step, sensitive, blot compatible.



## Order Today

Enjoy the benefits of SYPRO Orange. Order today or call to find out more about this unique stain. In the U.S., call 1-800-4BIORAD (1-800-424-6723) or contact your local Bio-Rad office.

Catalog No.	Product Description	Price
170-3120	<b>SYPRO Orange Concentrate,</b> 500 $\mu$ l, sufficient for 50–100 minigels	\$129.00
161-0332	<b>SDS-PAGE Standards for SYPRO Orange Staining,</b> broad range, 6.5–200 kd, 200 $\mu$ l (other size ranges available)	93.00

This product is offered for research purposes only. Coomassie is a registered trademark of ICI Organics, Inc. Polaroid is a registered trademark of Polaroid Corp. SYPRO is a registered trademark of, and the dye is manufactured by, Molecular Probes, Inc. Patents pending.



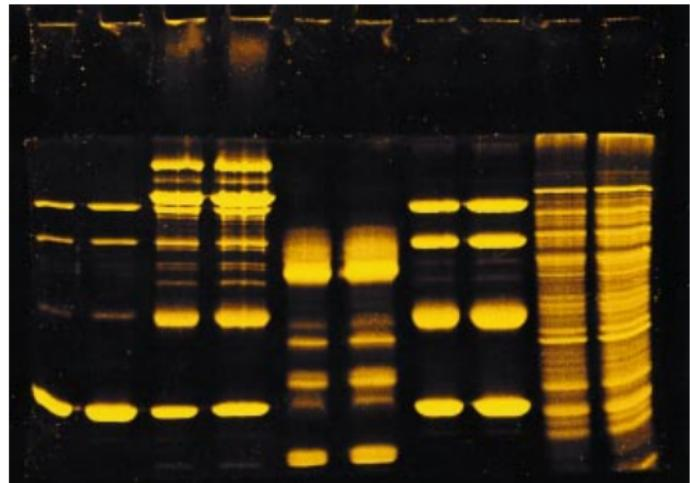
Bio-Rad  
Laboratories

Molecular  
Bioscience Group

U.S. (800) 4BIORAD California (510) 741-1000 Australia 02-9914-2800 Austria  
(1) 877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771 China (01) 2046622  
Denmark 39 17 9947 Finland 90 804 2200 France (1) 49 60 68 34 Germany 089  
31884-0 India 91-11-461-0103 Italy 02-21609 1 Japan 03-5811-6270 Hong Kong  
7893300 The Netherlands 0316-540661 New Zealand 09-443 3099 Singapore (65)  
272-9877 Spain (91) 661 70 85 Sweden 46 (0) 735 83 00 Switzerland 01-809 55 55  
United Kingdom 0800 181134

# New! SYPRO® Orange

## High-Sensitivity Fluorescent Protein Gel Stain



One 30-minute step, sensitive, blot compatible.



## Fast, Easy and Sensitive Staining

SYPRO Orange stain is a revolutionary breakthrough for staining proteins following gel electrophoresis. The SYPRO Orange reagent provides a reliable and reproducible fluorescent alternative to silver and Coomassie® staining. This unique dye is highly sensitive in the range of 1–10 ng per band when viewed photographically. SYPRO Orange requires only 30-60 minutes of staining in a single step.

Fluorescent protein bands are visualized by illuminating the gel with a 302 nm UV light box, similar to viewing ethidium bromide stained gels.

Photograph the fluorescent gel using Polaroid® black and white film at f5.6 for 5–15 sec, or record and analyze results using the Gel Doc™ 1000 system with 1–5 sec integration time.

## Compatible with Blotting

Unlike silver staining or Coomassie blue staining, stained gels are compatible with blotting or protein sequencing.

## Increase Productivity

SYPRO Orange stain increases productivity in SDS, native, and 2-D applications:

	SYPRO Orange	General Silver Stain	Coomassie Brilliant Blue
<b>Ease-of-use</b>	1 step	3–7 steps	2 steps
<b>Sensitivity</b>	1–10 ng/band	1–10 ng/band	30 ng/band
<b>Speed</b>	30–60 min	90 min	60 min
<b>Blot directly</b>	Yes	No	No

SYPRO Orange stain is virtually impossible to overdevelop and it will not stain DNA or RNA contaminants. It is also cost effective, because it is reusable.