
Criterion Staining and Blotting Trays

Catalog #
345-9920

For technical service, call your local Bio-Rad
office, or in the US, call 1-800-4BIORAD
(1-800-424-6723)

BIO-RAD

Introduction

Criterion™ staining/blotting trays are semi-disposable trays specifically designed for staining Criterion gels or performing western blot detection. Lightweight, compact, and stackable, the dimensions of the trays are optimized for Criterion gel staining and for optimal blot detection, providing a working volume of approximately 500 ml. Lids and spouts allow easy pouring without unnecessary gel handling.

Specifications

Material	PETG
Dimensions	15.8 cm x 9 cm (base) 17.3 cm x 11.1 cm (top) 3.6 cm (height)
Volume	500 ml

Stain Compatibility

BioSafe™ Coomassie stain
Coomassie Brilliant Blue, R-250
Coomassie Brilliant Blue, G-250
Bio-Rad silver stain
Bio-Rad Silver Stain Plus™ kit
SYPRO Ruby protein gel stain
SYPRO Orange stain
Copper
Zinc
Ethidium bromide

Staining and Blotting

For individual staining and blotting protocols, please refer to the corresponding stain or detection kit manual and proceed as indicated. After staining or blotting, discard remaining solutions by leaving tray lid on and tilting the tray so all liquid flows out over the pour spout. Trays may be reused multiple times; however, total number of uses depends on staining method.

Note: Microwaving or autoclaving of trays is not recommended.

Cleaning and Storage

Rinse trays thoroughly with deionized water and mild detergent. If necessary, wipe clean with ethanol for Coomassie staining or copper chloride for silver staining. Store stacked with or without lid.

This product is offered for research purposes only.

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

Bio-Rad Laboratories, Inc.
2000 Alfred Nobel Drive,
Hercules, CA 94547 USA
(510) 741-1000

1-800-424-6723

4110122 Rev A