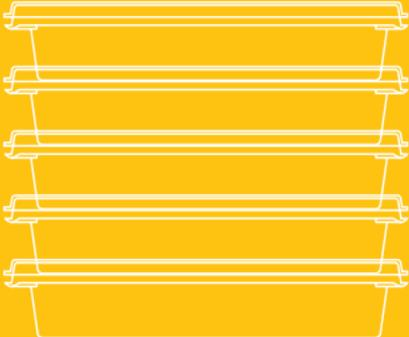


criterion.

A Tray for All Reasons

**Bio-Rad introduces convenient, semi-disposable,
stackable staining and blotting trays**

Tired of endless washing — but not ready to throw away your costly staining tray?
Worried about impurities left behind by previous staining? The new Criterion™
staining/blotting trays are low-cost, semi-disposable trays specifically designed
for staining one or two Criterion gels or performing western blot detection.
Leave tedious tray washing and oversized trays behind!



BIO-RAD

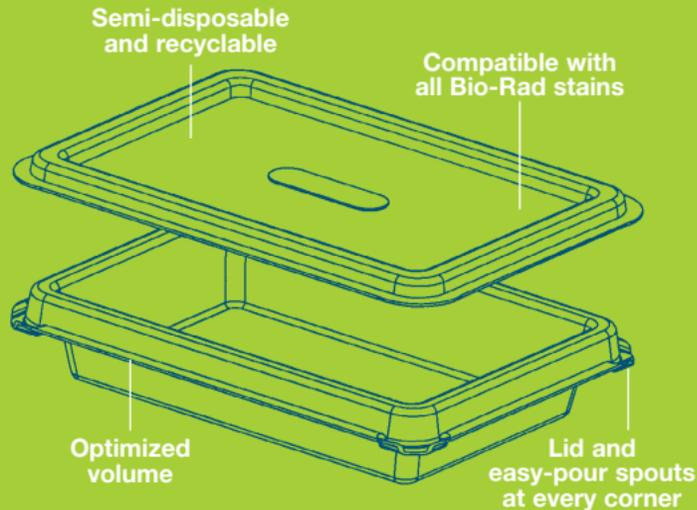
Stain and blot with the greatest of ease

These recyclable trays are lightweight, compact, and stackable. Their dimensions are optimized for Criterion gel staining and blot detection, providing a working volume of approximately 500 ml. Lids and spouts allow easy pouring without touching the gel.

Ordering Information

Catalog # Description

345-9920 Criterion Staining/Blotting Trays,
12 trays with lids



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 21 507 6191
Canada (905) 712-2771 **China** (86-21) 63052255 **Denmark** 45 44 52-1000 **Finland** 358 (0)9 804 2200 **France** 01 47 95 69 65 **Germany** 089 318 84-177
Hong Kong 852-2789-3300 **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** 52 5 534 2552 to 54 **The Netherlands** 0318-540666 **New Zealand** 64-9-4152280
Norway 47-23-38-41-30 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-2729877 **South Africa** 00 27 11 4428508 **Spain** 590 5200
Sweden 46 (0)8-55 51 27 00 **Switzerland** 061 717-9555 **United Kingdom** 0800-181134