

CRITERION™ BLOT WESTERN PROCESSOR KIT

CATALOG NUMBER 170-3985

Contents:

Outer tray and lid, 1
Criterion blot trays, 4
Tubing kits for use with one, two or four trays at a time
4 liter waste bottle with lid and fittings for aspiration

Application:

This kit adapts the Western Processor for the larger size Criterion precast gels, while still allowing you to develop one, two, or four blots at a time. A larger waste bottle is included in anticipation of larger volumes of reagents being used. For full operating instructions refer to the manual that came with your Western Processor (Bio-Rad part number 400-6145), and simply replace the tray and lid with the larger size tray and lid and use the tubing kits and waste bottle supplied with this kit.

Technical Assistance

For help, contact your local Bio-Rad representative. In the United States, call Technical Service at 1-800-4BIORAD.