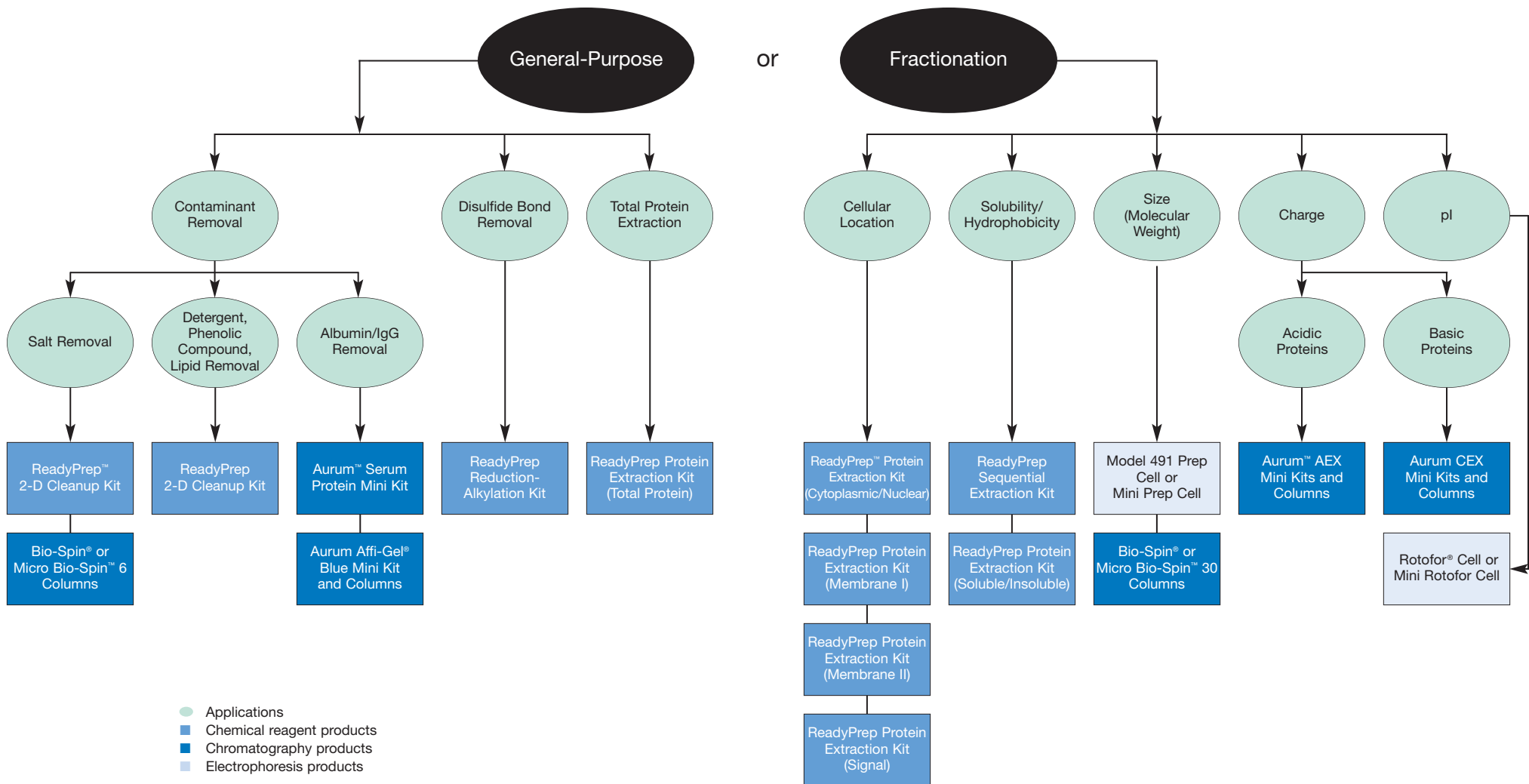


Product Selection Guide

Bio-Rad's comprehensive suite of 2-D sample preparation products.

These products may be used with virtually all sample types. They are all based on well-understood purification and fractionation principles and may be applied individually or in combination for effective sample preparation.



For more information, please visit us on the Web at www.expressionproteomics.com

Ordering Information

General-Purpose

Catalog # Description

Contaminant Removal: Salt Removal

163-2130 ReadyPrep 2-D Cleanup Kit
732-6227 Bio-Spin 6 Columns, in Tris buffer, 25
732-6228 Bio-Spin 6 Columns, in Tris buffer, 100
732-6002 Bio-Spin 6 Columns, in SSC buffer, 25
732-6213 Bio-Spin 6 Columns, in SSC buffer, 1,000
732-6221 Micro Bio-Spin 6 Columns, in Tris buffer, 25
732-6222 Micro Bio-Spin 6 Columns, in Tris buffer, 100
732-6225 Micro Bio-Spin 6 Columns, in Tris buffer, 1,000
732-6200 Micro Bio-Spin 6 Columns, in SSC buffer, 25
732-6201 Micro Bio-Spin 6 Columns, in SSC buffer, 100
732-6205 Micro Bio-Spin 6 Columns, in SSC buffer, 1,000

Contaminant Removal: Detergent, Phenolic Compound, and Lipid Removal

163-2130 ReadyPrep 2-D Cleanup Kit

Contaminant Removal: Albumin/IgG Removal

732-6713 Aurum Serum Protein Mini Kit, 2 pack
732-6701 Aurum Serum Protein Mini Kit, 10 pack
732-6712 Aurum Affi-Gel Blue Mini Kit, 2 pack
732-6708 Aurum Affi-Gel Blue Mini Columns, 25
732-6709 Aurum Affi-Gel Blue Mini Columns, 100

Disulfide Bond Removal

163-2090 ReadyPrep Reduction-Alkylation Kit

Total Protein Extraction

163-2086 ReadyPrep Protein Extraction Kit (Total Protein)

Fractionation

Catalog # Description

Cellular Location

163-2089 ReadyPrep Protein Extraction Kit (Cytoplasmic/Nuclear)
163-2088 ReadyPrep Protein Extraction Kit (Membrane I)
163-2084 ReadyPrep Protein Extraction Kit (Membrane II)
163-2087 ReadyPrep Protein Extraction Kit (Signal)

Solubility/Hydrophobicity

163-2100 ReadyPrep Sequential Extraction Kit
163-2085 ReadyPrep Protein Extraction Kit (Soluble/Insoluble)

Size (Molecular Weight)

170-2926 Model 491 Prep Cell and Pump, 100/120 V
170-2927 Model 491 Prep Cell and Pump, 220/240 V
170-2915 Mini Prep Cell With Reagent Starter Kit
170-2908 Mini Prep Cell Without Reagent Starter Kit
732-6231 Bio-Spin 30 Columns, in Tris buffer, 25
732-6232 Bio-Spin 30 Columns, in Tris buffer, 100
732-6006 Bio-Spin 30 Columns, in SSC buffer, 25
732-6223 Micro Bio-Spin 30 Columns, in Tris buffer, 25
732-6224 Micro Bio-Spin 30 Columns, in Tris buffer, 100
732-6226 Micro Bio-Spin 30 Columns, in Tris buffer, 1,000
732-6250 Micro Bio-Spin 30 Columns, RNase free, 25
732-6251 Micro Bio-Spin 30 Columns, RNase free, 100
732-6202 Micro Bio-Spin 30 Columns, in SSC buffer, 25
732-6203 Micro Bio-Spin 30 Columns, in SSC buffer, 100
732-6206 Micro Bio-Spin 30 Columns, in SSC buffer, 1,000

Charge: Acidic Proteins

732-6710 Aurum AEX Mini Kit, 2 pack
732-6705 Aurum AEX Mini Kit, 10 pack
732-6706 Aurum AEX Mini Columns, 25
732-6707 Aurum AEX Mini Columns, 100

Charge: Basic Proteins

732-6711 Aurum CEX Mini Kit, 2 pack
732-6702 Aurum CEX Mini Kit, 10 pack
732-6703 Aurum CEX Mini Columns, 25
732-6704 Aurum CEX Mini Columns, 100

Isoelectric Point

170-2986 Rotofor Purification System, 100/120 V
170-2987 Rotofor Purification System, 220/240 V
170-2988 Mini Rotofor Cell, 100/120 V, with starter kit
170-2989 Mini Rotofor Cell, 220/240 V, with starter kit