

Proteomics



Stellar Results

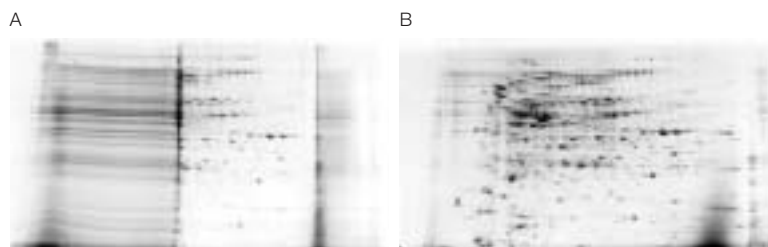
Get brilliant 2-D results with Bio-Rad's constellation of new ReadyPrep™ sample prep kits.



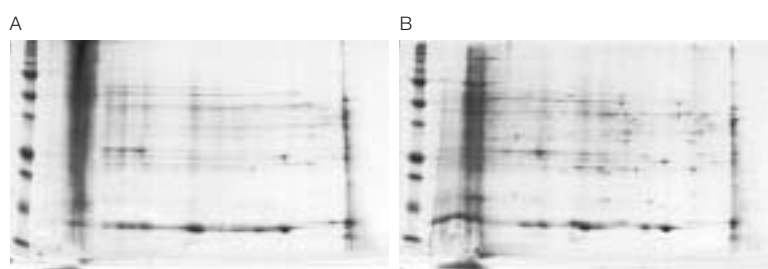
BIO-RAD

Get clear 2-D results.

One of the most critical steps in 2-D protein electrophoresis is sample preparation, which occurs even before the first-dimension separation begins. Many samples contain impurities that cloud the picture with streaking and spurious spots. Some samples require fractionation to enrich for low-abundance proteins. Bio-Rad's new ReadyPrep kits for sample cleanup and fractionation provide tested tools and proven protocols that put clear 2-D results within easy reach. For a closer look at ReadyPrep, go to www.bio-rad.com/ad/readyprep/



Effective detergent removal by the ReadyPrep 2-D cleanup kit leads to improved 2-D SDS-PAGE separation. *E. coli* extracts were spiked with 1% SDS. A, untreated sample; B, sample treated with the ReadyPrep kit.



Improved 2-D SDS-PAGE separation of mouse liver extract after treatment with the ReadyPrep reduction-alkylation kit. Both samples were applied by cup loading onto 11 cm ReadyStrip™ pH 7–10 IPG strips. A, reduced with 50 mM DTT in rehydration/sample buffer; B, treated with the ReadyPrep kit.

Application	ReadyPrep Kit	Extraction Method	# of Preps	Catalog #
General purpose	ReadyPrep 2-D cleanup kit	TCA-like precipitation	50	163-2130
	ReadyPrep reduction-alkylation kit	Reduction using TBP then alkylation with iodoacetamide	100	163-2090
	NEW ReadyPrep protein extraction kit (total protein)	Total protein solubilization	20	163-2086
Fractionation	ReadyPrep protein extraction kit (cytoplasmic/nuclear)	Organelle isolation followed by strongly chaotropic extraction buffer	50	163-2089
	ReadyPrep protein extraction kit (membrane I)	Temperature-dependent differential solubilization using Triton X-114	50	163-2088
	NEW ReadyPrep protein extraction kit (membrane II)	Carbonate buffer extraction	10	163-2084
	ReadyPrep protein extraction kit (signal)	Triton X-100 differential solubility	50	163-2087
	ReadyPrep sequential extraction kit	Differential solubility using SB 3-10	15	163-2100
	NEW ReadyPrep protein extraction kit (soluble/insoluble)	Differential solubility using ASB-14	20	163-2085

For more information about each of the kits, refer to bulletin 2934. Scientific poster reprints entitled Sample Preparation Solutions for Cellular Protein Fractionation are available by ordering bulletin 2961.

Triton is a trademark of Union Carbide.



**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 **Czech Republic** + 420 2 41 43 05 32 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **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** 55-52-00-05-20 **The Netherlands** 0318-540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** + 48 22 331 99 99 **Portugal** 351-21-472-7700 **Russia** 7 095 721 1404 **Singapore** 65-6415 3188 **South Africa** 00 27 11 4428508 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717-9555 **Taiwan** (8862) 2578-7189/2578-7241 **United Kingdom** 020 8328 2000