

Whole Gel Eluter

Bio-Rad Tech Note Summaries

Bulletin Title	#	Notes
Proteins		
Purification of Proteins From <i>Mycobacterium tuberculosis</i> by Simultaneous Electro-Elution on the Mini Whole Gel Eluter	2043	A procedure employing isoelectric focusing on a Rotofor cell, followed by SDS-PAGE and elution with the whole gel eluter
Purification of Inhibin Variants From Bovine Fluid Extract Using the Whole Gel Eluter and GS-670 Densitometer	2268	Native gel purification of affinity-enriched inhibin extract
Whole Gel Eluter Purification of a Functional Multiprotein DNA Replication Complex	2288	Native purification of a multienzyme complex. Elution is the final step in a purification scheme
Whole Gel Eluter Purification of Crosslinked High Density Lipoprotein and ApoA-1 and ApoA-2	2381	The whole gel eluter was used to recover HDL proteins. The resultant fractions were concentrated and passed over an AG11 column to remove free and bound SDS
Fractionation of <i>Bovicola ovis</i> Homogenates Using the Mini Whole Gel Eluter	2389	Two different elution buffer systems were compared for effectiveness of eluting a broad range of proteins
Combination of 2-D Gel and Liquid-Phase Electrophoretic Separations as Proteomic Tools in Neuroscience	2859	A summary of techniques incorporating the whole gel eluter, Model 491 prep cell, and Rotofor for prefractionation of complex protein mixtures in proteomics applications
Identification of Protein Vaccine Candidates From <i>Helicobacter pylori</i> Using a Preparative Two-Dimensional Electrophoretic Procedure and Mass Spectrometry	RP-0015	The Rotofor cell and whole gel eluter are used to separate detergent solubilized <i>H. pylori</i> membrane proteins prior to mass spectrometric analysis. This approach was successful in identifying a high percentage of membrane proteins that are typically underrepresented in 2-D PAGE gels
Proteomic study of a non-typable <i>Haemophilus influenzae</i>	RP-0025	Article describing a preparative 2-D separation involving first dimension separation on a Rotofor, followed by second dimension separation by SDS-PAGE and electroelution of fractions with the whole gel eluter. Use of this strategy in proteomic studies yields proteins not seen in traditional 2-D gel electrophoresis and in yields sufficient for characterization by MS
Nucleic Acids		
Isolating <i>Haloferax volcanii</i> Mature tRNA ^{Trp} -ProM From Total RNA Using the Whole Gel Eluter	2216	Mature tRNA is efficiently eluted from a polyacrylamide gel



**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 2527 3454
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 **Hungary** 36 1 455 8800 **India** (01-124)-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 12 700
Switzerland 061 717-9555 **Taiwan** (8862) 2578-7189/2578-7241 **United Kingdom** 020 8328 2000