

preparative electrophoresis

WHOLE GEL ELUTERS

SELECTED APPLICATIONS:

- EXPRESSED PROTEINS; QUICK RECOVERY OF FAT BANDS
- EPITOPE-TAGGED RECOMBINANT PROTEINS — SEPARATE FULL-LENGTH FUSION PROTEIN FROM PROTEOLYTIC FRAGMENTS CARRYING THE TAG
- PROTEIN DIGESTS — QUICKLY SEPARATE AND RECOVER FRAGMENTS
- RESOLVE THE DIGESTION, AND RECOVER THE FRAGMENTS
- mRNA, tRNA, 5S OR 7S RNA CAN BE EFFICIENTLY RECOVERED

Whole Gel Eluter and Mini Whole Gel Eluter — Purification as Simple as Electrophoresis.

Features of the Two Systems

The Whole Gel Eluter can recover any biomolecule that can be resolved on a slab gel. It should be useful in every lab that expresses recombinant proteins, or that uses electrophoresis to monitor a purification scheme. The mini Whole Gel Eluter is useful in any lab that uses mini gel apparatus, the Whole Gel Eluter can be placed in every lab that runs standard gels.

VERSATILE

- Elutes biomolecules from SDS- and native-PAGE, acid/urea gels, IEF acrylamide gels, agarose gels, slab gels up to 3 mm thick

QUICK

- 15 min–1 hr beyond running the resolving gel
- You can use the eluted protein by the time you Coomassie® stain a portion of the gel

HIGH RECOVERIES

- Recoveries of 70% or higher
- More efficient than electro-eluting an excised band
- A genuine advance over crush-and-soak method

PURITY

- Proteins resolved by more than 5 mm can be recovered in pure form
- Recombinant proteins can be quickly purified from proteolytic fragments



A Simple Concept for Purification

PROTEINS

Why not just pull that band right out of the gel? You run gels to monitor expression levels, and purification yields. Why not run an extra single lane gel and recover your protein in the time it takes to destain a Coomassie gel? The protein can be recovered in the buffer of your choice (choosing a buffer above the pI of the protein helps with the transfer). This method is ideal for inclusion body preps where your overexpressed protein can be dissolved in SDS, then eluted and refolded. This is also ideal where proteins can be resolved on native gels for and recovered as non-denatured proteins. Reaction mixtures can be quickly and easily separated. If your biomolecule will move through an electrical field, this quick and easy method will work for you!

NUCLEIC ACIDS

Nucleic acids separated on acid-urea PAGE, linear polymer acrylamide gels, PAGE, or agarose gels (less than 3 mm thick) can be successfully eluted with the whole gel eluter. If you can do an electrophoretic transfer to paper, you can transfer to liquid instead (transfer times may be 1 hr or greater).

BIO-RAD

How the Whole Gel Eluter works

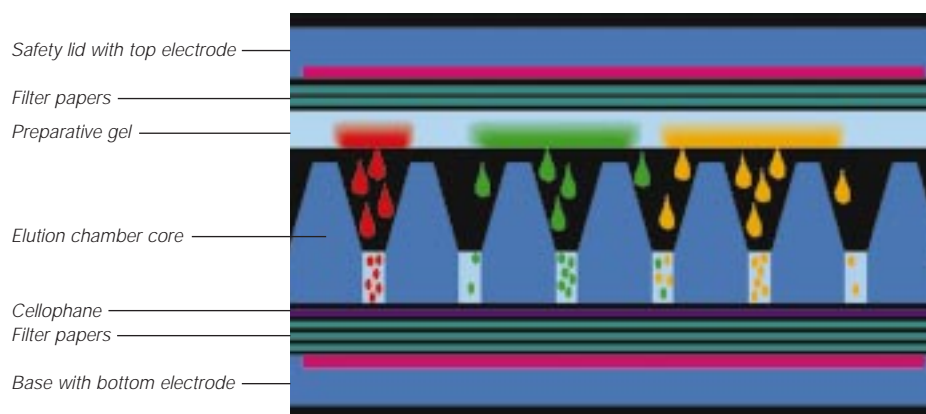
Place a whole preparative gel horizontally on the Whole Gel Eluter, preparative bands parallel to the channels of the eluter. Plate electrodes, which create a uniform electrical field, rapidly transfer RNA, DNA, or proteins from the gel into closely-spaced channels containing elution buffer. After the biomolecules are eluted into the corresponding channels, the fractions can be manually collected or immediately collected with the optional harvesting box attached to a vacuum source.

Specifications

	Whole Gel Eluter	Mini Whole Gel Eluter
FRACTIONS	30	14
GEL SIZE	14 x 16 cm or larger	6.5–5.5 cm or larger
MIN. GEL WIDTH	14 cm	5.5 cm
GEL THICKNESS	0.75–3.0 mm	0.75–3.0 mm
FRACTION VOL.	3.0 ml	0.5 ml
RUN TIME	15–20 min.	15–20 min.
POWER LIMIT	300 V/15 W	200 V/10 W
TOTAL BUFFER	1,000 ml	500 ml

Literature

Request the Whole Gel Eluter Technical Folder (Bulletin 1555C).



Proteins are electrophoretically eluted from the gel into precisely-spaced channels of the elution chamber core. Liquid fractions are then harvested.

Ordering information

Catalog #	Description
165-1256	Mini Whole Gel Eluter with Harvesting Box, includes Mini Whole Gel Eluter, Harvesting Box, all consumables for the first 25 uses, instructions
165-1255	Mini Whole Gel Eluter, includes Mini Whole Gel Eluter, all consumables for the first 25 uses, instructions
165-1261	Mini Harvesting Box
165-1287	Mini Whole Gel Eluter Consumables (complete set for 25 uses)
165-1251	Whole Gel Eluter with Harvesting Box, includes Whole Gel Eluter, Harvesting Box, all consumables for the first 25 uses, instructions
165-1250	Whole Gel Eluter, includes Whole Gel Eluter, all consumables for the first 25 uses, instructions
165-1260	Harvesting Box
165-1286	Whole Gel Eluter Consumables (complete set for 25 uses)

US Patent 08290993; Europe: 92905561.4; Japan: 504883/92; Canada: 2,130,751
Coomassie is a trademark of ICI Organics, Inc.

BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

Website www.bio-rad.com U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771
China 86-10-62051850 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France (01) 43 90 46 90 Germany 089 318 84-0
Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 Korea 82-2-3473-4460
The Netherlands 31 318-540666 New Zealand 64-9-4152280 Singapore 65-2729877 Spain (91) 661 70 85 Sweden 46 (0)8 627 50 00
Switzerland 01-809 55 55 United Kingdom 0800-181134