



290PR-02

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

EndotoxinOUT™

(Cat. # 786-366)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

EndotoxinOUT™ consists of 6% cross-linked agarose covalently linked to polymyxin B to bind and remove harmful pyrogens from a solution. Polymyxin B is a family, polymyxin B1 and B2, of antibiotics that bind to the negatively charged site of the lipid A portion of bacterial lipopolysaccharide layer neutralizing the endotoxic activity.

The covalent coupled agarose and polymyxin B is a stable matrix that resists leaching. An ideal product for the clean up of buffers, cell culture media, protein solutions, nucleic acid (DNA) samples and pharmacological components.

ITEM(S) SUPPLIED

Part #	Description	Size *
089E	EndotoxinOUT™	5 x 1mL
191E	EndotoxinOUT™ Regeneration Buffer	250mL
130W	Endotoxin Free Water	250mL

**EndotoxinOUT™ resin is supplied as a 50% slurry with 20% ethanol as a preservative.*

STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store refrigerated at 4°C, **DO NOT FREEZE**. This product is stable for 1 year at 4°C.

SPECIFICATIONS

- Capacity: ≥9995 endotoxin units (EU) removed by 1ml resin from 5ml test containing 10,000EU. ≥99.95% removal.
- Fractionation Range: 10-4,000kDa for proteins
- Bead Structure: 6% cross-linked agarose

IMPORTANT INFORMATION

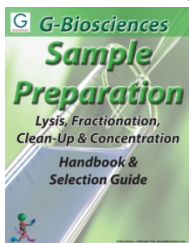
- Non specific binding of hydrophobic molecules may occur. To limit non-specific binding buffer all solutions to physiological pH. To inhibit ionic interactions, use a final concentration of 0.1-0.5M NaCl.
- Avoid the use of chaotropes and detergents as this interfere with binding to polymyxin sulfate.
- Some proteins, including BSA, bind tightly to endotoxins reducing their interaction and removal by the polymyxin sulfates. Increasing the volume of EndotoxinOUT™ resin to endotoxin containing sample may help.
- In some cases, the interaction between endotoxin and polymyxin sulfate is not inhibited, resulting in immobilization of the protein to the resin.
- Optimal performance of the EndotoxinOUT™ is best achieved under gravity flow chromatography, with an incubation period, allowing for more time for endotoxins to be in contact with polymyxin sulfate.

PROTOCOL

- Regenerate the EndotoxinOUT™ before and after every use.
 - Use endotoxin free solutions to prevent further endotoxin contamination.
 - Degas all solutions to prevent introducing air bubbles that inhibit column flow.
1. Allow the columns and reagents to equilibrate to room temperature.
 2. *Regeneration:* Wash the columns with 5ml EndotoxinOUT™ Regeneration Buffer, then wash with 5ml Endotoxin-Free Water.
 3. Equilibrate the column with 5ml of an appropriate endotoxin-free buffer or water.
 4. Apply the sample to the column and add the bottom then top cap. Incubate at room temperature for 30-60 minutes.
 5. Elute the sample with appropriate aliquots of endotoxin-free buffer or water. Repeat the elution 3-6 times and monitor elutions by absorbance at 280nm or an appropriate assay.
NOTE: Use extreme caution when handling samples to prevent contamination from poor handling, dirty glassware, etc.
 6. Repeat step 2 to regenerate the column and store in 25% ethanol at 4°C. The resin can be regenerated at least 10 times.

RELATED PRODUCTS

Download our Sample Preparation Handbook.



<http://info.gbiosciences.com/complete-protein-sample-preparation-handbook/>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 9/27/2016 TN



www.GBiosciences.com