



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

A Geno Technology, Inc. (USA) brand name

Gentle IgG Elution Buffer

(Cat. # 786-200, 786-201, 786-202)



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

INTRODUCTION

G-Biosciences Gentle IgG Elution Buffer is designed for the non-denaturing, high yield purification of antibodies from IgG affinity purification resins. The buffer is a near-neutral, high-salt conditions for elution, enabling even sensitive and labile antibodies (or other proteins) to be eluted from an affinity system without denaturation and inactivation. It is ideal for the purification of antibodies from Protein A, Protein G and Protein A/G resins.

ITEM(S) SUPPLIED

Cat. #	Description	Size
786-200	Gentle IgG Elution Buffer	100ml
786-201	Gentle IgG Elution Buffer	1L
786-202	Gentle IgG Elution Buffer	1gal

STORAGE CONDITION

The kit is shipped at ambient temperature. Upon arrival store the buffers at 4°C. Stable for 1 year when stored and used as recommended.

IMPOTANT INFORMAITON

Avoid using binding and wash buffers that contain phosphate ions (e.g., phosphate-buffered saline, PBS) because these will cause precipitation in the sample when the Gentle IgG Elution Buffer is applied. Use a non-phosphate binding and wash buffer that has the appropriate pH and ionic strength for the affinity interaction being used. Typical Ag/Ab binding interactions occur optimally at physiological conditions (pH 7.2-7.4, 150mM salt), and a solution with these parameters can be made with Tris, HEPES, MOPS or another non-phosphate buffer.

PROTOCOL: PURIFICATION OF IgG

Additional Item(s) Required

- Binding & Wash Buffer: Use a phosphate free buffer
- IgG Purification Resin
i.e. Immobilized Protein A, Protein G, Protein A/G or Protein L
- Neutralization Buffer: 1M Tris, pH8.0

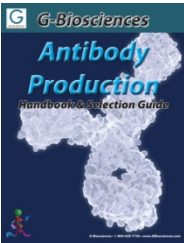
Procedure

1. Allow the buffers and resins to equilibrate to room temperature
2. Add an appropriate volume of Iresin to a suitable disposable column.
3. Equilibration Step: Wash the resin, by the addition of 5-10 column volumes (CV) of Binding/Wash Buffer. Allow wash/binding buffer to drain under gravity.

4. Gently apply the sample to the column by adding to the top of the resin. Do not disturb the gel bed.
Wash Step: Wash the column with 5-10CV of Binding/Wash Buffer or until the absorbance (280nm) of the flow through is near or at background levels.
5. Elution Step: Elute the immunoglobulins from the column by adding 5CV of elution buffer. Collect the eluate in 0.5-1ml fractions and immediately neutralize the elutions with 100µl Neutralization Buffer for every 1ml eluate.
6. Identify the immunoglobulin-containing fractions using a suitable protein assay or absorbance at 280nm.
7. Following elution, wash the resin with 5CV elution buffer, followed by at least 5CV of a suitable storage buffer containing a preservative. Store resin in storage buffer at 4°C.

RELATED PRODUCTS

Download our Antibody Production Handbook.



<http://info.gbiosciences.com/complete-Antibody-Production-handbook>

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

Last saved: 2/7/2018 CMH



www.GBiosciences.com