



572PR-01

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# IgG Binding, Wash & Elution Buffers

(Cat. # 786-203, 786-204, 786-205, 786-206, 786-544,  
786-545)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

G-Biosciences IgG Binding, Wash & Elution Buffers are designed for the non-denaturing, high yield purification of antibodies from IgG affinity purification resins. The buffers are ideal for the purification of antibodies from Protein A, Protein G and Protein A/G resins.

## ITEM(S) SUPPLIED

Cat. #	Description	Size
786-544	IgG Binding/ Wash Buffer	100ml
786-203	IgG Binding/ Wash Buffer	1L
786-204	IgG Binding/ Wash Buffer	1gal
786-545	IgG Elution Buffer	100ml
786-205	IgG Elution Buffer	1L
786-206	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.

## PROTOCOL: PURIFICATION OF IgG

### **Additional Item(s) Required**

- 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 IgG Purification Resin to a suitable disposable column.
3. Equilibration Step: Wash the resin, by the addition of 5-10 column volumes (CV) of IgG 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.

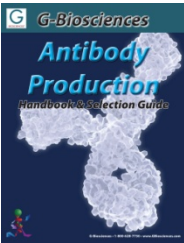
*NOTE: It is recommended to dilute serum samples at least 1:1 in the IgG Binding/Wash Buffer before applying to the resin. An alternative is to dialyze or perform a buffer exchange against IgG Binding/Wash Buffer.*

5. Wash Step: Wash the column with 5-10CV of IgG Binding/Wash Buffer or until the absorbance (280nm) of the flow through is near or at background levels.

6. **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.
7. Identify the immunoglobulin-containing fractions using a suitable protein assay or absorbance at 280nm.
8. 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](http://www.GBiosciences.com) or contact us.

Last saved: 2/7/2018 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)