



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

A Geno Technology, Inc. (USA) brand name

## DetergentOUT™ Tween® Gravity Flow Columns For Removal of Polysorbate (Tween®) Detergents

### KIT COMPONENTS

Cat. #	Description	Sample Size	Resin Volume	Size
786-2021	DetergentOUT™ Tween® Gravity flow column	6-10 ml	15 ml	10 columns
786-2022	DetergentOUT™ Tween® Gravity flow column	10-20 ml	30 ml	10 columns
786-2023	DetergentOUT™ Tween® Gravity flow column	50-100 ml	150 ml	1 column
786-2024	DetergentOUT™ Tween® Gravity flow column	50-100 ml	150 ml	5 columns



think proteins! think G-Biosciences!

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

## INTRODUCTION

Detergent-OUT™ Tween® resin has a high binding affinity for polysorbate detergents, commercially known as Tween® 20/80. The binding capacity is approximately 85mg polysorbate detergent per 1ml resin. The resin does not bind to protein allowing for a >95% protein recovery. DetergentOUT™ Tween® resin binds other commonly used detergents, but with lower affinity compared to DetergentOUT™ GBS10.

## STORAGE CONDITIONS

Shipped at ambient temperature. Store at 4°C, do not freeze.  
Storage buffer: 20% Methanol

## ADDITIONAL ITEMS NEEDED

- Collection tubes
- Equilibration Buffer

## PROCEDURE: DETERGENTOUT™ TWEEN® GRAVITY FLOW COLUMNS (15ML, 30 ML, 150 ML or LARGER COLUMNS)

### Column optimization:

Depending on sample type, it is recommended to optimize load capacity for each column. Load sample, elute small fraction size and analyze to establish load capacity. Generally, we recommend sample load not to exceed 50-60% of resin bed volume.

### Step by step procedure:

1. Remove the column from the Ziplock bag. Make sure the top and bottom caps are in a close position.
2. Open the top cap and remove stopper foams.
3. Place the column vertically with a clamp holder. Leave the column 30 minutes to equilibrate with the temperature.
4. Column contains 20% Methanol as a storage buffer. Drain it off completely.
5. Equilibrate the column with minimum 3 bead volume of the starting buffer when the pH and/or conductivity of the effluent is the same as the starting buffer.

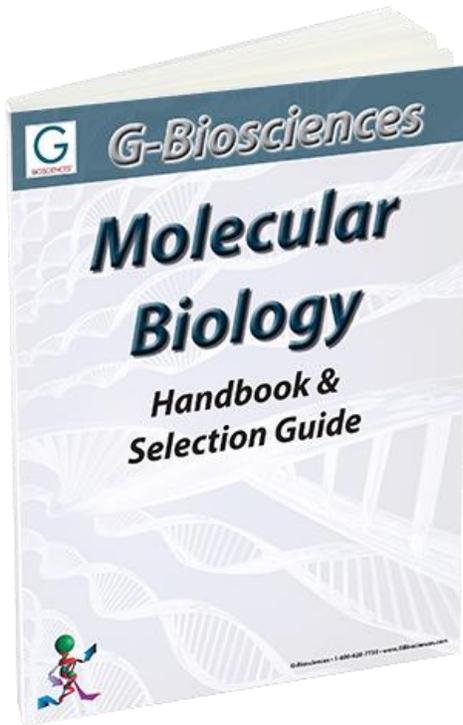
*NOTE: Do not equilibrate the columns with high concentration (>50mM) organic buffers, such as Tris buffer and other organic buffers as these may reduce the detergent binding capacity.*

6. Add the appropriate volume of detergent containing aqueous protein/peptide solution as indicated in the Kit Components. Remove the bottom cap and allow the sample to enter the resin.
7. Seal the column and incubate at room temperature for 2 minutes.
8. Elute and collect sample in clean collection tubes with an appropriate detergent free, aqueous buffer. Monitor elution with UV absorbance or an appropriate assay.

*Tween is a registered trademark of Uniqema, a business unit of ICI Americas, Inc.*

Download our Molecular Biology Handbook.

<http://info.gbiosciences.com/complete-molecular-biology-handbook/>



For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.