



**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®

## For Removal of Polysorbate (Tween®) Detergents

### KIT COMPONENTS

Cat. #	Description	Sample Size	Resin Volume	Size
786-2019	DetergentOUT™ Tween® (Maxi)	2-6 ml	10 ml	10 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®. 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.

## ADDITIONAL ITEMS NEEDED

- Collection tubes
- Equilibration Buffer (Deionized water)

## PREPARATION BEFORE USE

1. Prior to use, spin the DetergentOUT™ Tween® columns with the top and then bottom caps in place at 1,000xg for 1 minute to pellet the resin.
2. Mark one side of the column and ensure in all centrifugations the mark is facing outwards during centrifugation.

## PROCEDURE: DETERGENTOUT™ TWEEN® (MAXI) FOR 2-6 mL SAMPLE

1. Remove the bottom then top cap of the column. Place the column in a 50ml collection tube and centrifuge for 3 minutes at 1,000xg to remove the storage buffer.
2. Add 10ml Equilibration Buffer and then centrifuge for 2 minutes at 1,000xg. Discard the flow-through.

*NOTE: Detergent-OUT™ columns are supplied in deionized water. For optimal polysorbate detergent removal, equilibrate the column with deionized water. Columns may be equilibrated with other types of buffers, however, 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.*

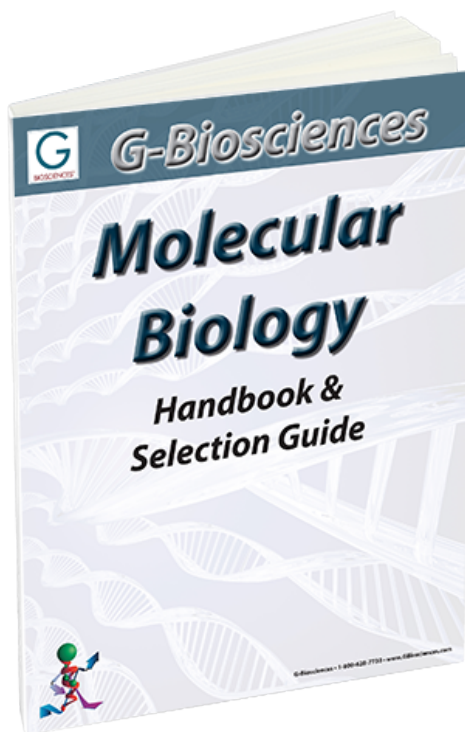
3. Repeat the equilibration step two more times.

4. Put on the bottom stopper and add 2-6 ml detergent containing aqueous protein/peptide solution. Put the top cap on and incubate at room temperature while shaking or inverting column (preferred), for 15 to 30 minutes.
5. Place column into a 50ml collection tube, centrifuge for 2 minutes at 1,500xg to collect the detergent-free sample.

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.