



## Certificate of Analysis / Data Sheet

**Product:** DetergentOUT Tween, Maxi

**Catalog No:** 786-2019

**Release Date:** 05/09/2023

**Size:** 10 columns

**Retest Date:** 04/29/2029

**Lot No:** 230105

**Storage:** 4 °C

### Specifications

**Appearance:** White resin

Test	Specification	Result
Binding Capacity	≥ 85mg Polysorbate detergent per ml resin	PASS
Protein Recovery	≥ 95%	PASS
Functional Test	Performs as per corresponding and/or companion protocol	PASS

This document has been produced electronically and is valid without signature

**Date of Print:** June 8, 2023

**G-Biosciences/Geno Technology, Inc.**

