



Certificate of Analysis / Data Sheet

Product: TCEP Reducing Resin

Catalog No: 786-822

Release Date: 7/27/20

Size: 1ml

Retest Date: 7/27/25

Lot No: 202707

Storage: 4°C

Specifications

CAS Number: N/A

Appearance: White, free of flocculent material

Test	Specification	Result
Reducing Capacity	≥ 8μmoles TCEP/ml of resin	20μmoles TCEP/ml resin

This document has been produced electronically and is valid without signature

Date of Print: August 18, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

