



Certificate of Analysis / Data Sheet

Product: TCEP-HCl

Catalog No: 786-874

Release Date: 02/03/2021

Size: 10g

Retest Date: 05/2023

Lot No: 210302

Storage: 4°C

Specifications

CAS Number: 51805-45-9

Appearance: White crystalline solid

Test	Specification	Result
Assay (NMR and HPLC)	> 98%	99.50%
Melting point	176.0 – 186.0°C	PASS
NMR (DMSO-d6)	Free of ammonium chloride	Conforms
Odor	Typical	Conforms
Solubility (water or DMSO)	Clear Solution	Conforms
Protease Activity	None Detected	Conforms

This document has been produced electronically and is valid without signature

Date of Print: February 19, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

