



Certificate of Analysis / Data Sheet

Product: TCEP-HCl

Catalog No: 786-874

Size: 10g

Lot No: 182012

Release Date: 01/04/2019

Retest Date: 04/20/2020

Storage: 4°C

Specifications

CAS Number: 51805-45-9

Appearance: White crystalline solid

Test	Specification	Result
Assay (NMR and HPLC)	> 98%	98.64%
Melting point	174 - 180°C	177.1 – 177.8°C
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: January 9, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

