



Certificate of Analysis / Data Sheet

Product: EDC (1-Ethyl-3-[3-dimethylaminopropyl]carbodiimide hydrochloride)
Catalog No: BC25-25 **Release Date:** 01/21/2019
Size: 25g **Retest Date:** 01/11/2020
Lot No: 191101 **Storage:** -20°C

Specifications

CAS Number: 25952-53-8

Appearance: White to off-white crystalline powder

Test	Specification	Result
Assay (Argentometric titration)	min 98%	99.9%
Melting Point (DSC)	110 - 114°C	112.2 to 112.4°C
Identity (IR spectrum)	Conforms to structure	Conforms
Clarity test	max 10NTU	2NTU
Water content (Karl Fischer)	max 0.5%	0.1%

This document has been produced electronically and is valid without signature

Date of Print: December 20, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

