



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** EDC; 1-ethyl-3-[3-dimethylaminopropyl]carbodiimide hydrochloride  
**Catalog #:** BC25-50  
**Quantity:** 50mg  
**Lot#:** 212801  
**MW:** 191.70  
**CAS#:** 25952-53-8

### Tests:

### Specification

Appearance	White to off white powder	Passed	Yes / <del>No</del>
Solubility	Faint yellow to yellow solution at 100mg/ml in water	Passed	Yes / <del>No</del>
Infrared Spectrum	consistent with structure	Passed	Yes / <del>No</del>
Carbon -	49.7 to 50.7	Passed	Yes / <del>No</del>
Nitrogen -	21.4 to 22.4	Passed	Yes / <del>No</del>

**Storage Temp:** Store at -20°C.

This document has been produced electronically and is valid without signature

**Date of Print:** February 11, 2021

### **G-Biosciences / Genotech**

9800 Page Avenue  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504

EDC C of A/Com, Rev 07.24.12/MM



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)