



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: 3,3',5,5'-Tetramethylbenzidine dihydrochloride (TMB HCl)

Catalog No: RC-309

Release Date: 10/5/21

Size: 100g

Retest Date: 10/5/23

Lot No: 210510

Storage: Desiccated at -20°C. Protect from light

Specifications

CAS Number: 207738-08-7

Appearance: White to off-white powder

Test	Specification	Result
Color of Solution (1% in methanol)	Colorless to light yellow/green	Light yellow
Turbidity (1% in methanol)	≤4.0NTU	0.2NTU
Purity (TLC)	≥97.0%	PASS

This document has been produced electronically and is valid without signature

Date of Print: October 13, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com