



Certificate of Analysis / Data Sheet

Product: 3,3',5,5'-Tetramethylbenzidine, dihydrochloride hydrate
Catalog No: RC-308 **Release Date:** 2/4/2019
Size: 10g **Retest Date:** 9/18/2019
Lot No: 191801 **Storage:** -20°

Specifications

CAS Number: 207738-08-7
Appearance: white powder

Test	Specification	Result
Color of Solution (1% in MeOH)	colorless to light yellow or light green-yellow	almost colorless
Turbidity (1% in MeOH)	≤ 4.0 NTU	0.3 NTU
Water Content (K. Fischer)	≤ 5.0%	0.3%
Sodium Hydroxide Titration (0.1 mol/L) (dry matter)	99.0 to 101.0%	9.7%
Sulfated Ash	≤ 0.1%	0.2%
FT-IR Spectrum	corresponds to reference spectrum	corresponds

This document has been produced electronically and is valid without signature

Date of Print: December 23, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

