



## Certificate of Analysis / Data Sheet

**Product:** Molecular Grade Water [Not DEPC Treated]

**Catalog No:** 786-292

**Release Date:** 10/29/2021

**Size:** 500ml

**Retest Date:** 10/29/2026

**Lot No:** 212910

**Storage:** Room Temp.

### Specifications

**CAS Number:** 7732-18-5

**Appearance:** Clear, colorless liquid

Test	Specification	Result
RNA Integrity Test (Overnight at 40°C)	PASS/FAIL	PASS
DNase / RNase / Protease Activity	No Activity	PASS
Sterility	No Growth (37°C Overnight)	PASS
Conductivity	< 1.0ppm, < 1.5µS	PASS
pH	4.50-5.00	PASS

This document has been produced electronically and is valid without signature

**Date of Print:** December 21, 2021

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP

