



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

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

**Catalog No:** 786-1774

**Release Date:** 12/06/2021

**Size:** 55 Gallons

**Retest Date:** 12/06/2026

**Lot No:** 210612

**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:** January 13, 2022

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

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)