



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Molecular Grade Water [Not DEPC Treated]

Catalog No: 786-293

Release Date: 10/11/2021

Size: 1L

Retest Date: 10/11/2026

Lot No: 211110

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 10, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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