



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

**Product:** Endotoxin Free Water (<0.0050 EU/ml)

**Catalog #:** 786-670

**Size:** 500ml

**Lot #:** 212910

**Re-test date:** 10/2026

**Appearance:** Clear Liquid

### QC Tests:

Endotoxin level <0.0050 EU/ml	Passed	Yes / <del>Ne</del>
RNA Integrity Test (overnight at 40°C)	Passed	Yes / <del>Ne</del>
RNase / DNase Activity - None Detected	Passed	Yes / <del>Ne</del>
Protease Activity – None Detected	Passed	Yes / <del>Ne</del>
Sterility Test (overnight at 37°C)	Passed	Yes / <del>Ne</del>
Conductivity (<1.0ppm, <1.5 µS,)	Passed	Yes / <del>Ne</del>
pH range (4.5- 5.0)	Passed	Yes/ <del>Ne</del>

**Storage:** Store at Room Temp. It can also be stored at 4°C for long term storage.

*This document has been produced electronically and is valid without signature*

*For Research Use Only; Not for Cell Culture or Human Use*

**Date of Print:** November 12, 2021

### **G-Biosciences / Genotech**

9800 Page Avenue

St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504

Endotoxin Free Water 786-671, C of A/Com#3/Rev 07.17.12/MM



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)