



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:** HPLC Grade Water

**Catalog #:** 786-1798

**Size:** 4L

**Lot #:** 200511

**Re-test date:** 11/2025

**Appearance:** Clear Liquid

### Test(s):

RNase / DNase / Protease Activity - None Detected	Passed	Yes / <del>No</del>
Sterility Test (overnight at 37°C)	Passed	Yes / <del>No</del>
Conductivity (<1.0ppm, <1.5 µS)	Passed	Yes / <del>No</del>
pH range (~5.0- 6.0 ± 0.25)	Passed	Yes/ <del>No</del>
Absorbance (1.3mAU @ 254nm)	Passed	Yes/ <del>No</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 4, 2021

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

9800 Page Avenue  
St. Louis, MO 63132, USA  
Ph: 314-991-6034, Fax: 314-991-1504



think proteins! think G-Biosciences!

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