



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

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

Catalog #: 786-671

Size: 1L

Lot #: 212809

Appearance: Clear Liquid

Re-test date: 09/2031

QC Test (s):

Endotoxin level <0.0050 EU/ml	Passed	Yes / No
RNA Integrity Test (overnight at 40°C)	Passed	Yes / No
RNase / DNase Activity - None Detected	Passed	Yes / No
Protease Activity – None Detected	Passed	Yes / No
Sterility Test (overnight at 37°C)	Passed	Yes / No
Conductivity (<1.0ppm, <1.5 µS,)	Passed	Yes / No
pH range (4.5- 5.0)	Passed	Yes/ No

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: October 8, 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#3/07.17.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com