



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

Certificate of Analysis

Product: Ultrapure Elite™ Water

Catalog #: 786-1789

Size: 1L

Lot #: 212106

Appearance: Clear Liquid

Re-test date: 06/2031

QC Test (s):

Endotoxin level <0.0050 EU/ml	Passed	Yes / No
RNase / DNase / Protease Activity - None Detected	Passed	Yes / No
Sterility Test (overnight at 37°C)	Passed	Yes / No
Conductivity (<1.0ppm, <1.5 µS _c)	Passed	Yes / No
pH range (4.5- 5.0)	Passed	Yes/ No
Meets All ACS Grade Specifications	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: 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