



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

Certificate of Analysis

Product: Molecular Grade Water
[Not DEPC Treated]

Catalog #: 786-292

Size: 500ml

Lot #: 191607

Appearance: Clear Liquid

Test(s):

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 (~5.0- 6.0 ± 0.25)	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: September 20, 2019

Expiry Date: July 16, 2024

G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

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

Molecular Grade Water 786-292, C of A/Com#3



think proteins! think G-Biosciences!

www.GBiosciences.com



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

Rev 03.16.15/IA



think proteins! think G-Biosciences!

www.GBiosciences.com