



Certificate of Analysis

Product: [20X] SSPE

Catalog #: R022

Size: 1 Gallon

Lot #: 202307

Re-test date: 07/2025

Tests:

Appearance - Clear Liquid	Passed	Yes / No
Specific gravity test	Passed	Yes / No
pH test	Passed	Yes / No

Storage: Store at room temperature

For Research Use Only

This document has been produced electronically and is valid without signature

Date of Print: August 7, 2020

G-Biosciences / Genotech

9800 Page Avenue

St. Louis, MO 63132, USA

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

www.GBiosciences.com

20X SSPE Cat# R022 C of A/Com#3

Rev 01.17.13/MM



think proteins! think G-Biosciences!

www.GBiosciences.com