



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

Certificate of Analysis

Product: [20X] SSC

Catalog #: 786-023

Lot #: 171805

Re-test date: 05/2022

Size: 500ml

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

20X SSC Cat# 786-023 C of A/Com#3
Rev 03.24.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com