



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Sodium Chloride [5M]

Catalog No: R006

Release Date: 7/30/20

Size: 100ml

Retest Date: 7/30/25

Lot No: 203007

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: Clear, colorless liquid

Test	Specification	Result
pH @ RT	7.5 – 9.0	8.93

This document has been produced electronically and is valid without signature

Date of Print: January 14, 2022

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com