



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Sodium Chloride [5M]

**Catalog No:** R006

**Release Date:** 12/03/20

**Size:** 100ml

**Retest Date:** 12/03/25

**Lot No:** 200312

**Storage:** Room Temp.

### Specifications

**CAS Number:** N/A

**Appearance:** Clear, colorless liquid

Test	Specification	Result
pH @ RT	6.9 – 8.5	8.03

This document has been produced electronically and is valid without signature

**Date of Print:** December 21, 2020

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)