



**G-Biosciences** ♦

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

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Hybridization Neutralizing Solution

**Catalog No:** R018

**Release Date:** 11/2/20

**Size:** 1 gallon

**Retest Date:** 11/2/24

**Lot No:** 200211

**Storage:** Room Temp.

### Specifications

**CAS Number:** N/A

**Appearance:** Clear, colorless solution

Test	Specification	Result
pH	7.35 – 7.45	7.41

This document has been produced electronically and is valid without signature

**Date of Print:** November 6, 2020

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

Rev 10/25/18 BP



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