



**G-Biosciences** ♦

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

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** TMAC Hybridization Solution [3X], pH 8.0  
**Catalog No:** 786-538 **Release Date:** 01/25/21  
**Size:** 1L **Retest Date:** 01/21/25  
**Lot No:** 212501 **Storage:** Room Temp.

### Specifications

**CAS Number:** N/A

**Appearance:** Clear, colorless solution. Heat to 37°C if precipitated.

Test	Specification	Result
pH	7.90 – 8.10	8.06

This document has been produced electronically and is valid without signature

**Date of Print:** February 19, 2021

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

Rev 10/25/18 BP



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