



**G-Biosciences** ♦

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

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Yeast tRNA (10mg/ml)

**Catalog No:** 786-058

**Release Date:** 8/6/20

**Size:** 0.5ml

**Retest Date:** 8/6/22

**Lot No:** 200608

**Storage:** -20°C

### Specifications

**CAS Number:** N/A

**Appearance:** Clear liquid

Test	Specification	Result
Integrity of tRNA	Meets specs.	PASS
RNase Test	None detected	PASS

This document has been produced electronically and is valid without signature

**Date of Print:** August 18, 2020

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

Rev 10/25/18 BP



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