



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: DNAOUT™

Catalog No: 786-75

Release Date: 01/28/21

Size: 1L

Retest Date: 01/28/27

Lot No: 212801

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: Clear, colorless liquid

Test	Specification	Result
pH	> 12.00	12.94

This document has been produced electronically and is valid without signature

Date of Print: February 11, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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