



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: SDS-PAGE Gel Fixing Solution

Catalog No: 786-236

Release Date: 11/5/19

Size: 1L

Retest Date: 11/5/24

Lot No: 190511

Storage: Room Temp.

Specifications

Appearance: Clear solution

Test	Specification	Result
pH	2.24-3.24	2.69
Gel fixing test	Pass test	PASS

This document has been produced electronically and is valid without signature

Date of Print: November 18, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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