



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: HIS Elution Buffer, pH 8.0

Catalog No: 786-543

Release Date: 11/21/19

Size: 100ml

Retest Date: 11/21/22

Lot No: 192111

Storage: Room Temp.

Specifications

Appearance: Clear, colorless liquid

Test	Specification	Result
pH	7.95 – 8.05	8.00

This document has been produced electronically and is valid without signature

Date of Print: December 2, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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