



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Protein G IgG Binding Buffer

Catalog No: 786-1601

Release Date: 7/21/20

Size: 1L

Retest Date: 7/21/23

Lot No: 202107

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: Clear, colorless liquid

Test	Specification	Result
pH	4.9 – 5.1	5.05

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