



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Antigen Retrieval Solution, Citrate (pH 6.0) [50X]

Catalog No: 786-1260

Release Date: 03/17/2023

Size: 500ml

Retest Date: 03/15/2027

Lot No: 231603

Storage: 4°C

Specifications

CAS Number: Not applicable

Appearance: Clear, colorless to light yellow liquid

Test	Specification	Result
pH	5.8 – 6.2	PASS

This document has been produced electronically and is valid without signature

Date of Print: June 2, 2023

G-Biosciences/Geno Technology, Inc.

Rev 02/13/2022 DB



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