



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: RIPA Lysis and Extraction Buffer

Catalog No: 786-723

Release Date: 11/25/19

Size: 1L

Retest Date: 11/25/22

Lot No: 192511

Storage: Room Temp.

Specifications

Appearance: Clear, Colorless liquid

Test	Specification	Result
pH	6.5-8.5	7.62

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