



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Nonidet® P-40 Substitute

Catalog No: 786-512

Release Date: 8/12/2021

Size: 1 Liter

Retest Date: 12/24/2022

Lot No: 211008

Storage: Room Temp.

Specifications

CAS Number: 9016-45-9

Appearance: Clear, colorless

Test	Specification	Result
Appearance	Clear, colorless	Passed
Solubility (5% in H ₂ O)	Clear, colorless	Passed

This document has been produced electronically and is valid without signature

Date of Print: January 28, 2022

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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