



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Nonidet P-40

Catalog No: 786-512

Release Date: 9/22/20

Size: 1L

Retest Date: 9/22/26

Lot No: 202209

Storage: Room Temp.

Specifications

CAS Number: 9016-45-9

Appearance: Colorless to slightly yellow, viscous liquid

Test	Specification	Result
Density (20C)	1.05-1.07 g/ml	1.06 g/ml
Refractive Index	1.48-1.52	1.49
Solubility (5% in water)	PASS/FAIL	PASS

This document has been produced electronically and is valid without signature

Date of Print: September 29, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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