



Certificate of Analysis

Product: Nonidet[®] P-40 Substitute [Nonylphenyl-polyethylene glycol]

Catalog #: DG501C

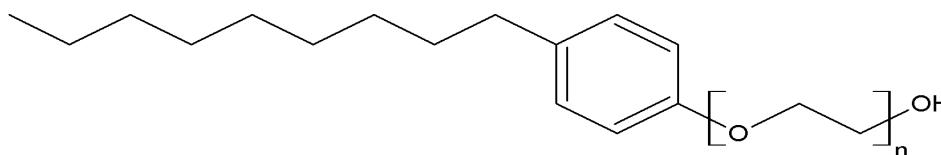
Size : 1 x 50ml

Lot#: 200911

Re-test date: 11/2025

Detergent Type: Non-ionic detergent

Structure:



Mol. Formula: C₁₅H₂₄O [C₂H₄O]_n **Mol Weight:** 573 (for n=8)

Form: 10% aqueous solution (w/v) **Absorbance** (254nm): 0.14 (0.05% w/v)

Typical Analysis:

Aldehyde content < 200µM

Peroxide content (as H₂O₂) < 50µM

Application: A commonly used non-ionic detergent for solubilizing membrane proteins during isolation of membrane-protein complexes. Ultra low aldehyde and peroxide concentrations reduces the effects of peroxidase and carbonyl compounds that negatively interact with membrane proteins.

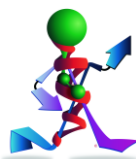
Critical micelle concentration (CMC): Approx 0.05-0.3mM (25°C)

Storage: Store at 4°C, protected from light. It is sterile and packaged under inert gas and stable for 1 year as supplied.

This document has been produced electronically and is valid without signature

Date of Print: November 20, 2020

Nonidet P-40 Cat# DG501 C of A/Com#3/12.14.09/Rev IA



think proteins! think G-Biosciences!

www.GBiosciences.com