



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: 2D-Detergent Nonidet[®] P-40 Substitute [Nonylphenyl-polyethylene glycol]
10% Aqueous Solution, Purified, Sealed vials under argon gas

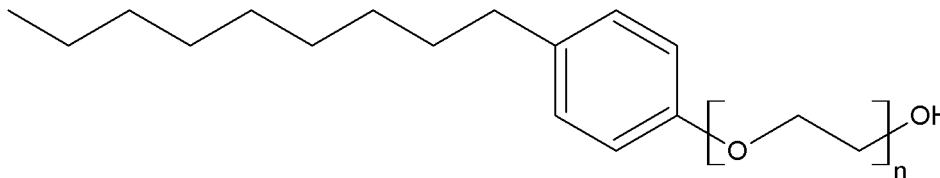
Catalog #: DG902

Size: 10 x 10ml

Lot#: 101202

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 < 100μM

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.

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

This document has been produced electronically and is valid without signature

Date of Print: October 31, 2011

2D Detergent Nonidet P-40 Cat# DG902 C of A/Com#3
Rev 10.31.11/IA



think proteins! think G-Biosciences!

www.GBiosciences.com