



## Certificate of Analysis

**Product:** Nonidet<sup>®</sup> P-40 Substitute [Nonylphenyl-polyethylene glycol]

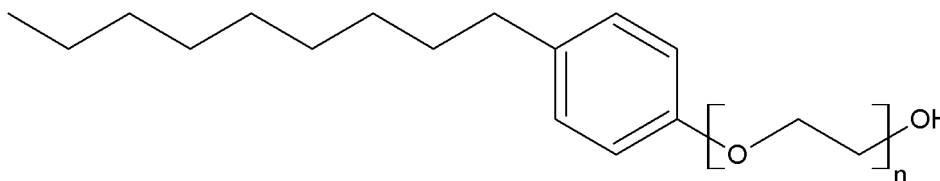
**Catalog #:** DG514C

**Size :** 1 x 100ml

**Lot#:** 200911

**Detergent Type:** Non-ionic detergent

**Structure:**



**Mol. Formula:**  $C_{15}H_{24}O [C_2H_4O]_n$

**Mol Weight:** 573 (for n=8)

**Form:** 10% aqueous solution (w/v)

**Absorbance (254nm):** 0.14 (0.05% w/v)

**Typical Analysis:**

**Aldehyde content** < 200 $\mu$ M

**Peroxide content (as H<sub>2</sub>O<sub>2</sub>)** < 50 $\mu$ M

**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.

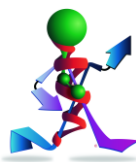
**Critical micelle concentration (CMC):** approx 0.05-0.3mM (25°C)

**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# DG514 C of A/Com#3  
Rev 06.13.11/IA



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)