



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

Certificate of Analysis

Product: Triton[®] X-100 [Octylphenolpoly(ethyleneglycolether)_x]

Catalog #: DG007; DG008

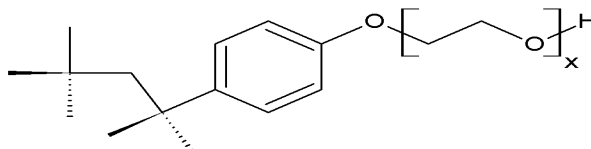
Size: 5 x 10ml; 10 x 10ml

Lot#: 200110

Re-test date: 10/2025

Detergent Type: Non-ionic detergent

Structure:



Mol. Formula: C₃₄H₆₂O₁₁ for x =10

Mol Weight: 647 (for x=10)

Form: 10% aqueous solution (w/v)

Absorbance (254nm): 0.16 (0.05% w/v)

Typical Analysis:

Aldehyde content < 100μM

Peroxide content (as H₂O₂) < 50μM

Application: It is one of the most commonly used non-ionic detergents 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 0.2 x 10⁻³M (25°C)

Cloud Point: 65°C

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

This document has been produced electronically and is valid without signature

Date of Print: October 20, 2020

**Triton is a registered trademark of Union Carbide Corp.*

Triton-X100, DG008 C of A/Com#3
Rev 07.02.12/IA



think proteins! think G-Biosciences!

www.GBiosciences.com