



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

Certificate of Analysis

Product: 2-D Detergent™ Triton® X-100, 10% Aqueous Solution
[Purified and Sealed vials under Argon gas]

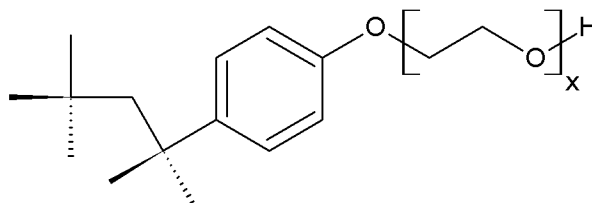
Catalog #: DG907

Size: 5 x 10ml

Lot#: 180811

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.

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

This document has been produced electronically and is valid without signature

Date of Print: February 21, 2019

2D Detergent Triton X-100 Cat# DG907 C of A/Com#3
Rev 09.08.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com