



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

CAS#: 9002-93-1

Detergent Type: Non-ionic detergent

Mol. Formula: C₃₄H₆₂O₁₁ for x = 10

Mol Weight: 647 (for x=10)

Form: 10% aqueous solution (w/v)

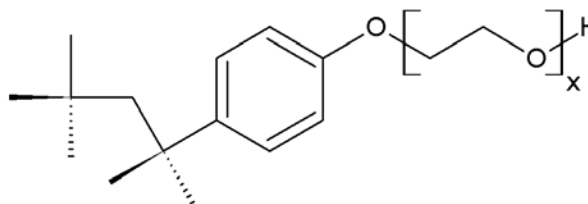
Absorbance (254nm): 0.16 (0.05% w/v)

Critical micelle concentration (CMC): Approx. 0.2 x 10⁻³M (25°C)

Cloud Point: 65°C

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.

Storage: Store at 2-8°C, protected from light. It is Sterile, packaged under inert gas and stable for 1 year.



<u>Test</u>	<u>Specification</u>	<u>Result</u>
Conductivity	≤15.0μS)	14.0 μS
Aldehyde concentration	≤100μM	0 μM
Peroxide concentration	≤50μM	0 μM
Detergent concentration	9.0-11%	10%
Sterility		Passed

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

This document has been produced electronically and is valid without signature

Date of Print: February 28, 2019
Rev 2/28/2019 CMH

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



think proteins! think G-Biosciences!

www.GBiosciences.com