



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## Certificate of Analysis

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

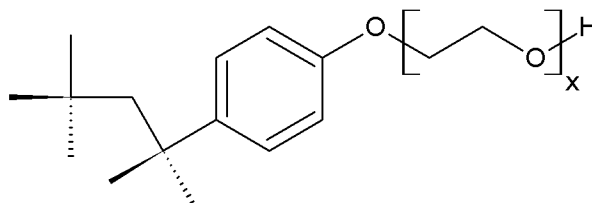
**Catalog #:** DG907

**Size:** 5 x 10ml

**Lot#:** 102211

**Detergent Type:** Non-ionic detergent

**Structure:**



**Mol. Formula:** C<sub>34</sub>H<sub>62</sub>O<sub>11</sub> 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 H<sub>2</sub>O<sub>2</sub>) < 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<sup>-3</sup>M (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 9, 2011

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