

Product Information



TBS-Polysorbate Stock Solution (10X, pH 7.4)

Item No. 600222

Contents: This bottle contains 250 mM Tris-base, pH 7.4, 1.37 M sodium chloride, containing 27 mM KCl, 0.4 mM phenol red, and 1.0% (v/v) polysorbate.

Stability: ≥ 1 year at 4°C

Laboratory Procedures

This bottle contains a concentrated solution of filtered TBS-polysorbate. It is ready to use as supplied. For long term storage, we suggest that this product be stored at 4°C; it should be stable for at least one year.

To prepare for use, dilute the concentrated TBS-polysorbate to 10 L by adding 1 part stock with 9 parts UltraPure water (*i.e.*, 1 L stock into 9L UltraPure water). If the concentrated TBS-polysorbate has been frozen, it may remain slightly gelatinous or contain crystals even after thawing. This is normal and will disappear upon dilution. The bottle should be rinsed with the diluted TBS-polysorbate to ensure that all the salts are removed. Diluted TBS-polysorbate may be stored for several weeks at room temperature; the pH of this solution will be ≈ 7.4 .

CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/600222

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/24/2014