

PRODUCT INFORMATION

(PEO)3-monoamine

Item No. 30426

CAS Registry No.: 6338-55-2
Formal Name: 2-[2-(2-aminoethoxy)ethoxy]-ethanol

Synonyms: NSC 40760, T₃EGMA, Triethylene glycol monoamine, Triglycolamine

MF: C₆H₁₅NO₃

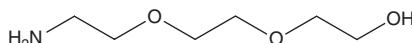
FW: 149.2

Purity: ≥98%

Supplied as: A liquid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(PEO)3-monoamine is supplied as a liquid. A stock solution may be made by dissolving the (PEO)3-monoamine in the solvent of choice, which should be purged with an inert gas. (PEO)3-monoamine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (PEO)3-monoamine in these solvents is approximately 15, 10, and 30 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of (PEO)3-monoamine can be prepared by directly dissolving the liquid in aqueous buffers. The solubility of (PEO)3-monoamine in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(PEO)3-monoamine is a poly(ethylene oxide) (PEO) blocking reagent that prevents nonspecific adsorption of analytes in applications using covalent protein coupling such as ELISA.¹ It reduces the adsorption of IgG when coated on carboxylic-terminated surfaces.

Reference

1. Frederix, F., Bonroy, K., Reekmans, G., *et al.* Reduced nonspecific adsorption on covalently immobilized protein surfaces using poly(ethylene oxide) containing blocking agents. *J. Biochem. Biophys. Methods* **58**(1), 67-74 (2004).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/25/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM