

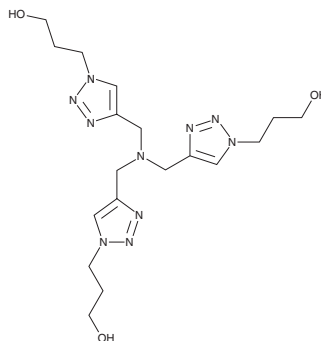
PRODUCT INFORMATION



THPTA

Item No. 35670

CAS Registry No.: 760952-88-3
Formal Name: 4,4',4''-[nitrilotris(methylene)]tris-1H-1,2,3-triazole-1-propanol
MF: $C_{18}H_{30}N_{10}O_3$
FW: 434.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

THPTA is supplied as a solid. A stock solution may be made by dissolving the THPTA in the solvent of choice, which should be purged with an inert gas. THPTA is soluble in DMSO.

A stock solution may be made by dissolving the THPTA in water. We do not recommend storing the aqueous solution for more than one day.

Description

THPTA is a ligand for copper(I)-catalyzed azide/alkyne cycloaddition (CuAAC).¹ It protects biomolecules from hydrolysis by copper(II) byproducts and against oxidation by various copper, oxygen, or ascorbate-derived radicals and peroxides.

Reference

1. Presolski, S.I., Hong, V.P., and Finn, M.G. Copper-catalyzed azide-alkyne click chemistry for bioconjugation. *Curr. Protoc. Chem. Biol.* **3**(4), 153-162 (2011).

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, 03/04/2024

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