

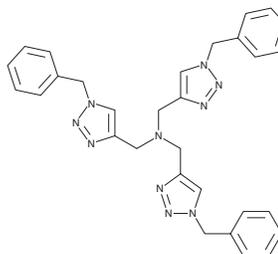
# PRODUCT INFORMATION



## TBTA

Item No. 18816

**CAS Registry No.:** 510758-28-8  
**Formal Name:** tris[(1-benzyl-1H-1,2,3-triazol-4-yl)methyl]amine  
**Synonyms:** Click Tag™ TBTA,  
Tris(benzyltriazolylmethyl)amine  
**MF:** C<sub>30</sub>H<sub>30</sub>N<sub>10</sub>  
**FW:** 530.6  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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

TBTA is supplied as a crystalline solid. A stock solution may be made by dissolving the TBTA in the solvent of choice. TBTA is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of TBTA in these solvents is approximately 30 mg/ml.

TBTA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TBTA should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TBTA has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

TBTA is a tertiary amine with three 1,2,3-triazole groups. It complexes with, and stabilizes, copper(I) to accelerate azide-alkyne cycloadditions, as used in click chemistry.<sup>1,2</sup>

### References

1. Hong, V., Udit, A.K., Evans, R.A., *et al.* Electrochemically protected copper(I)-catalyzed azide-alkyne cycloaddition. *ChemBioChem* **9(9)**, 1481-1486 (2008).
2. Kuang, G.-H., Michaels, H.A., Simmons, J.T., *et al.* Chelation-assisted, copper(II)-acetate-accelerated azide-alkyne cycloaddition. *J. Org. Chem.* **75(19)**, 6540-6548 (2010).

#### 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, 08/24/2023

#### 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