

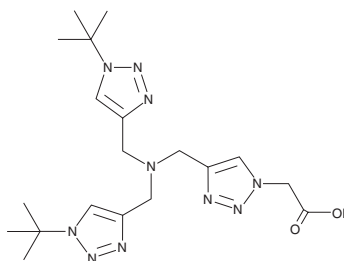
PRODUCT INFORMATION



BTAA

Item No. 41089

CAS Registry No.: 1334179-85-9
Formal Name: 4-[[bis[[1-(1,1-dimethylethyl)-1H-1,2,3-triazol-4-yl]methyl]amino]methyl]-1H-1,2,3-triazole-1-acetic acid
MF: C₁₉H₃₀N₁₀O₂
FW: 430.5
Purity: ≥98%
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

BTAA is supplied as a solid. A stock solution may be made by dissolving the BTAA in the solvent of choice, which should be purged with an inert gas. BTAA is soluble (10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

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 BTAA can be prepared by directly dissolving the solid in aqueous buffers. BTAA is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

BTAA is a click chemistry copper (I) (Cu(I))-stabilizing ligand and catalyst.¹ It displays an increased rate of Cu(I)-catalyzed azide-alkyne cycloaddition (CuAAC) in a propargyl alcohol and 3-azido-7-hydroxycoumarin reaction, as well as an increased yield, compared to the click chemistry ligands BTES, THPTA (Item No. 35670), and TBTA (Item No. 18816) in a fluorogenic assay. BTAA has been used with biotin-azide in CuAAC for the labeling and detection of azide-tagged glycoconjugates in Jurkat cells, zebrafish, and bacteria.²

References

1. Besanceney-Webler, C., Jiang, H., Zheng, T., *et al.* Increasing the efficacy of bioorthogonal click reactions for bioconjugation: a comparative study. *Comparative Study* **50(35)**, 8051-8056 (2011).
2. Besanceney-Webler, C., Jiang, H., Wang, W., *et al.* Metabolic labeling of fucosylated glycoproteins in Bacteroidales species. *Bioorg. Med. Chem. Lett.* **21(17)**, 4989-4992 (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, 07/02 /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