

PRODUCT INFORMATION



Methyltetrazine-PEG₅-NHS ester

Item No. 44734

CAS Registry No.: 1802907-92-1

Formal Name: 3-[2-[2-[2-[2-[4-(6-methyl-1,2,4,5-tetrazin-3-yl)phenoxy]ethoxy]ethoxy]ethoxy]ethoxy]propanoic acid, 2,5-dioxo-1-pyrrolidinyl ester

Synonym: MTzS

MF: C₂₄H₃₁N₅O₉

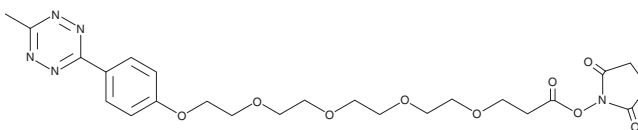
FW: 533.5

Purity: ≥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

Methyltetrazine-PEG₅-NHS ester is supplied as a solid. A stock solution may be made by dissolving the methyltetrazine-PEG₅-NHS ester in the solvent of choice, which should be purged with an inert gas. Methyltetrazine-PEG₅-NHS ester is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Methyltetrazine-PEG₅-NHS ester is a click chemistry building block.¹ It has been used to induce protein-DNA cross-linking and as a starting material in the synthesis of clickable antibodies.^{1,2}

References

1. Hayes, O.G., Partridge, B.E., and Mirkin, C.A. Encoding hierarchical assembly pathways of proteins with DNA. *Proc. Natl. Acad. Sci. U.S.A.* **118(40)**, e2106808118 (2021).
2. Komatsu, T., Kyo, E., Ishii, H., *et al.* Antibody clicking as a strategy to modify antibody functionalities on the surface of targeted cells. *J. Am. Chem. Soc.* **142(37)**, 15644-15648 (2020).

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, 10/27/2025

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