

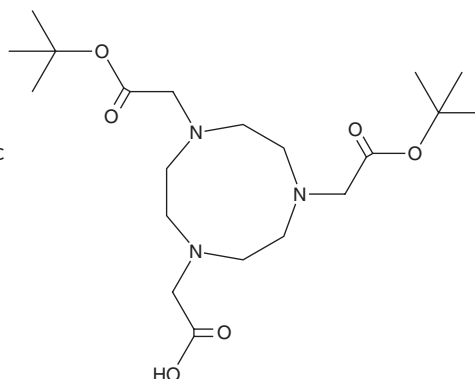
PRODUCT INFORMATION



NOTA-*bis*(*t*Bu ester)

Item No. 42964

CAS Registry No.: 1161415-28-6
Formal Name: hexahydro-1H-1,4,7-triazonine-1,4,7-triacetic acid, 1,4-*bis*(1,1-dimethylethyl) ester
Synonym: N(*t*EB)₂
MF: C₂₀H₃₇N₃O₆
FW: 415.5
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

NOTA-*bis*(*t*Bu ester) is supplied as a solid. A stock solution may be made by dissolving the NOTA-*bis*(*t*Bu ester) in the solvent of choice, which should be purged with an inert gas. NOTA-*bis*(*t*Bu ester) is soluble (≥10 mg/ml) in ethanol and 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 NOTA-*bis*(*t*Bu ester) can be prepared by directly dissolving the solid in aqueous buffers. NOTA-*bis*(*t*Bu ester) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

NOTA-*bis*(*t*Bu ester) is a derivative of the macrocyclic chelator NOTA.¹ It has been used in the synthesis of albumin-dimerizing conjugates that prevent peptide proteolysis *in vivo* and HER2-specific compounds and bombesin receptor 2-specific peptides chelated to radioactive metals for tumor imaging *in vivo*.¹⁻³

References

1. Tian, R., Zhu, S., Zeng, Q., *et al.* An albumin sandwich enhances *in vivo* circulation and stability of metabolically labile peptides. *Bioconjug. Chem.* **30**(6), 1711-1723 (2019).
2. Tolmachev, V., Altai, M., Sandström, M., *et al.* Evaluation of a maleimido derivative of NOTA for site-specific labeling of affibody molecules. *Bioconjug. Chem.* **22**(5), 894-902 (2011).
3. Varasteh, Z., Velikyan, I., Lindeberg, G., *et al.* Synthesis and characterization of a high-affinity NOTA-conjugated bombesin antagonist for GRPR-targeted tumor imaging. *Bioconjug. Chem.* **24**(7), 1144-1153 (2013).

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, 04/04/2025

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