

PRODUCT INFORMATION



tris-NTA (trifluoroacetate salt)

Item No. 28789

Formal Name: ($\alpha^1S, \alpha^4S, \alpha^8S$)-11-(6-amino-1-oxohexyl)- $\alpha^1, \alpha^4, \alpha^8$ -tris[bis(carboxymethyl)amino]- $\delta^1, \delta^4, \delta^8$ -trioxo-1,4,8,11-tetraazacyclotetradecane-1,4,8-tripentanoic acid, trifluoroacetate salt

Synonym: tris-Nitrilotriacetic Acid

MF: $C_{43}H_{72}N_8O_{22} \cdot XCF_3COO$

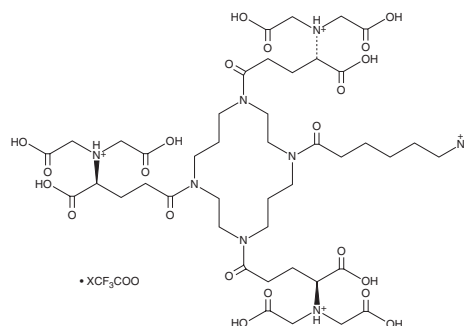
FW: 1,053.1

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

tris-NTA (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the tris-NTA (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

tris-NTA is a His-tagged protein ligand.^{1,2} It has been used alone, or conjugated to fluorophores, for the purification and detection of His-tagged proteins.

References

- Huang, Z., Hwang, P., Watson, D.S., et al. Tris-nitrilotriacetic acids of subnanomolar affinity toward hexahistidine tagged molecules. *Bioconjug. Chem.* **20(8)**, 1667-1672 (2009).
- Glembockyte, V., Wieneke, R., Gatterdam, K., et al. Tris-N-nitrilotriacetic acid fluorophore as a self-healing dye for single-molecule fluorescence imaging. *J. Am. Chem. Soc.* **140(35)**, 11006-11012 (2018).

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, 09/18/2024

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