

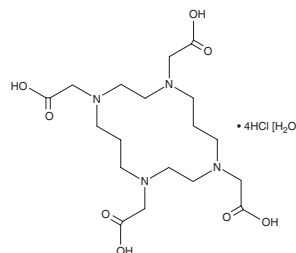
PRODUCT INFORMATION



TETA (hydrochloride hydrate)

Item No. 43750

CAS Registry No.: 314041-07-1
Formal Name: 1,4,8,11-tetraazacyclotetradecane-1,4,8,11-tetraacetic acid, tetrahydrochloride hydrate
MF: $C_{18}H_{32}N_4O_8 \cdot 4HCl [H_2O]$
FW: 596.3
Purity: $\geq 95\%$
Supplied as: A crystalline 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

TETA (hydrochloride hydrate) is supplied as crystalline solid. Aqueous solutions of TETA (hydrochloride hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. TETA (hydrochloride hydrate) is soluble (≥ 10 mg/ml) (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

TETA is a macrocyclic chelator.¹ It has been conjugated to radioligands for use in PET imaging of neuroendocrine tumors and for tissue distribution studies *in vivo*.

Reference

1. Lewis, J.S., Srinivasan, A., Schmidt, M.A., *et al.* *In vitro* and *in vivo* evaluation of ^{64}Cu -TETA-Tyr³-octreotate. A new somatostatin analog with improved target tissue uptake. *Nucl. Med. Biol.* **26**(3), 267-273 (1999).

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, 08/13/2025

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