

PRODUCT INFORMATION



CAY10799

Item No. 39619

CAS Registry No.: 1423622-48-3
Formal Name: N-[[[(1S)-2-(1,1-dimethylethoxy)-1-[[4-hydroxyphenyl)methyl]-2-oxoethyl]amino]carbonyl]-L-glutamic acid 1,5-bis(1,1-dimethylethyl) ester
Synonym: Di-tert-butyl (((S)-1-(tert-butoxy)-3-(4-hydroxyphenyl)-1-oxopropan-2-yl)carbamoyl)-L-glutamate

MF: C₂₇H₄₂N₂O₈

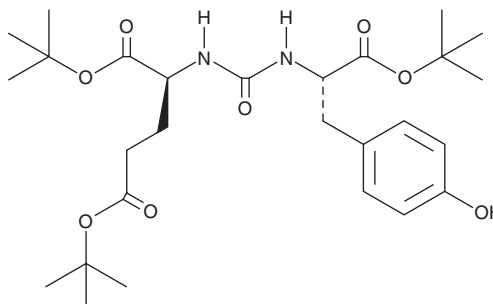
FW: 522.6

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

CAY10799 is supplied as a solid. A stock solution may be made by dissolving the CAY10799 in the solvent of choice, which should be purged with an inert gas. CAY10799 is soluble in methanol, acetonitrile, and chloroform.

Description

CAY10799 is a precursor in the synthesis of prostate-specific membrane antigen (PSMA) ligands and fluorescent probes that have been used in the detection of PSMA in LNCaP and PC3 cells.¹

Reference

1. Sengupta, S., Krishnan, M.A., Pandit, A., *et al.* Tyrosine-based asymmetric urea ligand for prostate carcinoma: Tuning biological efficacy through *in silico* studies. *Bioorg. Chem.* **91**, 103154 (2019).

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, 12/18/2023

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