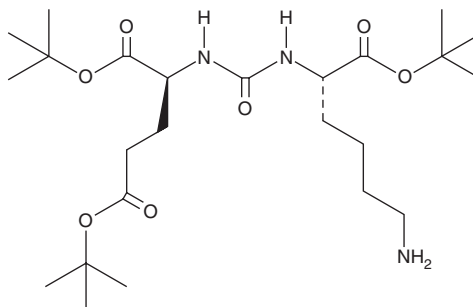


PRODUCT INFORMATION

Di-*tert*-butyl (((S)-6-amino-1-(*tert*-butoxy)-1-oxohexan-2-yl)carbamoyl)-L-glutamate

Item No. 39616

CAS Registry No.: 1025796-31-9
Formal Name: N-[[[(1S)-5-amino-1-[(1,1-dimethylethoxy)carbonyl]pentyl]amino]carbonyl]-L-glutamic acid, 1,5-bis(1,1-dimethylethyl) ester
MF: C₂₄H₄₅N₃O₇
FW: 487.6
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

Di-*tert*-butyl (((S)-6-amino-1-(*tert*-butoxy)-1-oxohexan-2-yl)carbamoyl)-L-glutamate is supplied as a solid. A stock solution may be made by dissolving the di-*tert*-butyl (((S)-6-amino-1-(*tert*-butoxy)-1-oxohexan-2-yl)carbamoyl)-L-glutamate in water. We do not recommend storing the aqueous solution for more than one day.

Description

Di-*tert*-butyl (((S)-6-amino-1-(*tert*-butoxy)-1-oxohexan-2-yl)carbamoyl)-L-glutamate is a synthetic intermediate that has been used in the synthesis of prostate-specific membrane antigen-targeted radioconjugates with anticancer activity.¹

Reference

1. Santos, J.F., Braz, M.T., Raposinho, P., *et al.* Synthesis and preclinical evaluation of PSMA-targeted ¹¹¹In-radioconjugates containing a mitochondria-tropic triphenylphosphonium carrier. *Mol. Pharm.* **21**(1), 216-233 (2024).

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, 01/23/2024

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