

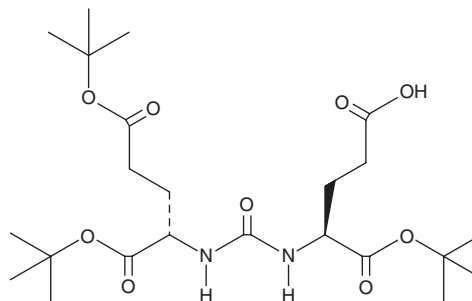
PRODUCT INFORMATION



DUPA(OtBu)-OH

Item No. 39617

CAS Registry No.: 1026987-94-9
Formal Name: N-[[[(1S)-3-carboxy-1-[(1,1-dimethylethoxy)carbonyl]propyl]amino]carbonyl]-L-glutamic acid, 1,5-bis(1,1-dimethylethyl) ester
Synonym: DUPA(OtBu)₃-OH
MF: C₂₃H₄₀N₂O₉
FW: 488.6
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

DUPA(OtBu)-OH is supplied as a solid. A stock solution may be made by dissolving the DUPA(OtBu)-OH in the solvent of choice, which should be purged with an inert gas. DUPA(OtBu)-OH is soluble in acetonitrile.

Description

DUPA(OtBu)-OH is a precursor in the synthesis of conjugates of the prostate-specific membrane antigen (PSMA) ligand 2-[3-(1,3-dicarboxypropyl)ureido]pentanedioic acid (DUPA).^{1,2} It has been used in the synthesis of PSMA-targeted radio- and optical imaging agents and prodrugs with anticancer activity.

References

1. Kularatne, S.A., Wang, K., Santhapuram, H.-K.R., *et al.* Prostate-specific membrane antigen targeted imaging and therapy of prostate cancer using a PSMA inhibitor as a homing ligand. *Mol. Pharm.* **6(3)**, 780-789 (2009)
2. Kularatne, S.A., Venkatesh, C., Santhapuram, H.-K.R., *et al.* Synthesis and biological analysis of prostate-specific membrane antigen-targeted anticancer prodrugs. *J. Med. Chem.* **53(21)**, 7767-7777 (2010).

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, 05/02/2024

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