

PRODUCT INFORMATION



MACTIDE Azide

Item No. 43997

Contents: This vial contains 1 mg of lyophilized peptide trifluoroacetic acid salt
Formal Name: (6S,9S,12R,15S,18R,23R,26S,28aS,33aS)-18-(2-(6-((S)-3-(4-(2λ⁴-triazol-1,2-dien-1-yl)phenyl)-2-aminopropanamido)hexanamido)-acetamido)-12-(4-aminobutyl)-6,26-di((S)-sec-butyl)-N-((S)-1-((S)-2-(((S)-1-(((S)-1,6-diamino-1-oxohexan-2-yl)amino)-1-oxopropan-2-yl)amino)-2-oxoethyl)carbamoyl)pyrrolidin-1-yl)-3-hydroxy-1-oxopropan-2-yl)-15-((R)-1-hydroxyethyl)-9-(hydroxymethyl)-5,8,11,14,17,25,28,33-octaoxo-octacosahydro-1H,5H,22H-dipyrrolo[2,1-j:2',1'-m][1,2]dithia[5,8,11,14,17,20,23,26]octaazacyclononacosine-23-carboxamide

Synonyms: UNO-related Cyclic Peptide,
CD206 Cyclic Peptide Ligand Azide

Peptide Sequence: 4-Azido-Phe-(Ahx)-G[CTKSIPPIC]-SPGAK

MF: C₇₇H₁₂₆N₂₃O₂₀S₂

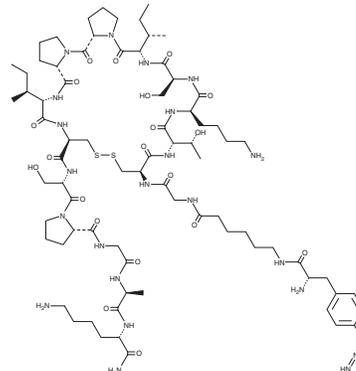
FW: 1,757.9

Form: A solid

Storage: -20°C (as supplied)

Stability: ≥3 years

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



Laboratory Procedures

MACTIDE azide is soluble in water, buffer, and ethanol up to 20 mg/ml. Used for the delivery of DBCO-labeled molecules to CD206. We do not recommend storing the aqueous solution for more than one day.

Description

MACTIDE azide is a clickable form of MACTIDE, a cyclic peptide that binds CD206 on tumor-associated macrophages (TAMs).¹

Reference

1. Lepland, A., Peranzoni, E., Haljasorg, U., *et al.* Peptide-drug conjugate for therapeutic reprogramming of tumor-associated macrophages in breast cancer. *Adv. Sci. (Weinh)* **12**(10), e2410288 (2025).

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, 09/04/2025

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