

# PRODUCT INFORMATION



## Azide-UNO Peptide

Item No. 43998

**Contents:** This vial contains 1 mg of lyophilized peptide trifluoroacetic acid salt.

**Formal Name:** (S)-2-((S)-2-((S)-2-((S)-2-((S)-2-((S)-1-((6-((S)-3-(4-(2(4-triaza-1,2-dien-1-yl)phenyl)-2-aminopropanamido)hexanoyl)-L-seryl)pyrrolidine-2-carboxamido)acetamido)propanamido)-6-aminohexanamido)-3-methylbutanamido)-5-guanidinopentanimidic acid

**Synonym:** CD206 Peptide Ligand Azide

**Peptide Sequence:** 4-Azido-Phe-Ahx-SPGAKVR

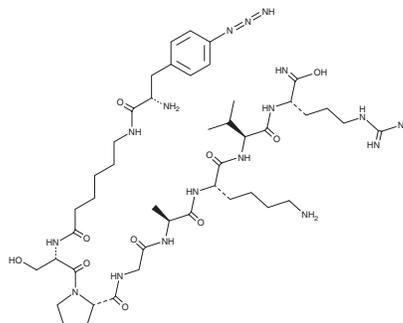
**MF:** C<sub>45</sub>H<sub>76</sub>N<sub>17</sub>O<sub>10</sub>

**FW:** 1,015.2

**Form:** A solid

**Storage:** -20°C

**Stability:** ≥3 years



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

### Laboratory Procedures

Azide-UNO peptide 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

Azide-UNO peptide is a clickable form of UNO peptide, a peptide that binds CD206 on tumor-associated macrophages (TAMs).<sup>1</sup>

### Reference

1. Leland, 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