

PRODUCT INFORMATION



LDV-K(azide) (trifluoroacetate salt)

Item No. 45188

Formal Name: (S)-4-(((S)-1-(((S)-1-amino-6-azido-1-oxohexan-2-yl)amino)-3-methyl-1-oxobutan-2-yl)amino)-3-((S)-4-methyl-2-(2-(4-(3-(*o*-tolyl)ureido)phenyl)acetamido)pentanamido)-4-oxobutanoic acid, trifluoroacetate salt

Synonyms: 2-(4-(3-(*o*-tolyl)ureido)phenyl)acetyl-Leu-Asp-Val-Lys(N₃), PUPA-LDV azide, PUPA-Leu-Asp-Val-Lys(azide)

Peptide Sequence: PUPA-LDV-K(azide) (where PUPA = phenylureidophenylacetyl)

MF: C₃₇H₅₂N₁₀O₈ • XCF₃COOH

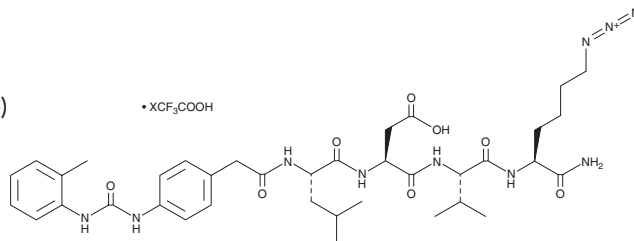
FW: 764.9

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

LDV-K(azide) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the LDV-K(azide) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. LDV-K(azide) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

LDV-K(azide) is a functionalized peptide containing an LDV motif for binding integrin α4β1 and a modified version of LDV-azide.¹ LDV-K(azide) can be conjugated to alkyne-containing molecules via click chemistry.

Reference

1. Komoriya, A., Green, L.J., Mervic, M., *et al.* The minimal essential sequence for a major cell type-specific adhesion site (CS1) within the alternatively spliced type III connecting segment domain of fibronectin is leucine-aspartic acid-valine. *J. Biol. Chem.* **266**(23), 15075-15079 (1991).

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, 05/27/2026

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