

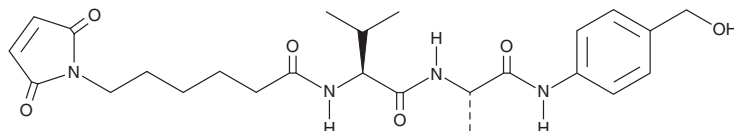
PRODUCT INFORMATION



Mc-Val-Ala-PAB

Item No. 43201

CAS Registry No.: 1870916-87-2
Formal Name: N-[6-(2,5-dihydro-2,5-dioxo-1H-pyrrol-1-yl)-1-oxohexyl]-L-valyl-N-[4-(hydroxymethyl)phenyl]-L-alaninamide
MF: $C_{25}H_{34}N_4O_6$
FW: 486.6
Purity: $\geq 98\%$
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

Mc-Val-Ala-PAB is supplied as a solid. A stock solution may be made by dissolving the Mc-Val-Ala-PAB in the solvent of choice, which should be purged with an inert gas. Mc-Val-Ala-PAB is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Mc-Val-Ala-PAB is a peptide linker molecule that has been used in the synthesis of antibody-drug conjugates (ADCs).^{1,2} It contains a maleimidocaproyl (Mc) group that can be conjugated to an antibody and a p-aminobenzyl (PAB) spacer that allows the peptide to be linked to active compounds, such as anticancer agents or antibiotics. Mc-Val-Ala-PAB cleavage *in vitro* and *in vivo* by cathepsin B allows the active compound to be released at the target site. It has been used in the synthesis of ADCs containing the antimitotic compound monomethyl auristatin E (MMAE; Item No. 16267) or the antibiotic KRM-1657.

References

1. Li, D., Sun, X., Li, Y., *et al.* AGCM-22, a novel cetuximab-based EGFR-targeting antibody-drug-conjugate with highly selective anti-glioblastoma efficacy. *Bioorg. Med. Chem.* **102**, 117657 (2024).
2. Fan, S., Bai, Y., Li, Q., *et al.* Novel antibody-antibiotic conjugate using KRM-1657 as payload eliminates intracellular MRSA *in vitro* and *in vivo*. *Bioorg. Chem.* **150**, 107532 (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, 04/09/2025

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