

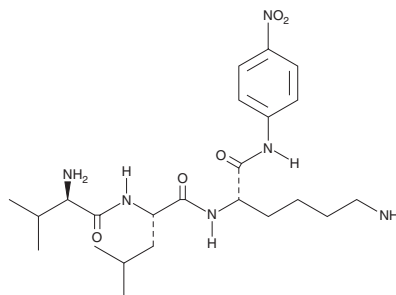
PRODUCT INFORMATION



D-Val-Leu-Lys-pNA

Item No. 36580

CAS Registry No.: 62354-43-2
Formal Name: D-valyl-L-leucyl-N-(4-nitrophenyl)-L-lysineamide
Synonyms: D-Val-Leu-Lys-*p*-nitroanilide,
D-VLK-*p*NA, S 2251
MF: C₂₃H₃₈N₆O₅
FW: 478.6
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 315 nm
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

D-Val-Leu-Lys-*p*NA is supplied as a solid. A stock solution may be made by dissolving the D-Val-Leu-Lys-*p*NA in water. We do not recommend storing the aqueous solution for more than one day.

Description

D-Val-Leu-Lys-*p*NA is a colorimetric substrate for plasmin.^{1,2} Plasmin binds and hydrolyzes D-Val-Leu-Lys-*p*NA to release *p*-nitroanilide (*p*NA), which can be quantified by colorimetric detection at 405 nm as a measure of plasmin activity.

References

1. Gohda, K., Fujimori, K., Teno, N., *et al.* Synthetic substrates specific to activated plasmin can monitor the enzymatic functional status *in situ* in breast cancer cells. *Chem. Biol. Drug Des.* **83**(1), 52-57 (2014).
2. Leite, J.A.S., Montoya, C.A., Loveday, S.M., *et al.* Heat-treatments affect protease activities and peptide profiles of ruminants' milk. *Front. Nutr.* **8**, 626475 (2021).

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, 09/18/2024

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