

Ac-VEID-pNA

Caspase substrate

Colorimetric substrate for caspase-6 ($K_m=30\mu\text{M}$; $K_{cat}/K_m=168'000\text{ M}^{-1}\text{sec}^{-1}$) and related cysteine proteases. Sequence is based on the lamin A site of cleavage. Cleavage of the substrate by a caspase releases pNA, which can be detected by its absorbance at 405nm ($\epsilon=10,500\text{ M}^{-1}\text{cm}^{-1}$).

Ordering Information

[Order Online »](#)

BML-P437-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Caspase-6 substrate (chromogenic)

Appearance White to off-white powder.

Formulation Lyophilized.

Identity Determined by MS.

MW 636.6

Purity ≥95% (HPLC)

Sequence N-acetyl-Val-Glu-Ile-Asp-pNA (p-nitroanilide)

Solubility Soluble at 1 mg/ml in water with 5% ammonium hydroxide.

Source Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)