

Ac-WEHD-pNA

Substrate for caspase-1, -4, -5 and -14

Colorimetric caspase substrate. Combinatorial studies have found the WEHD sequence to be optimal for cleavage of tetrapeptide substrate by caspase-1, caspase-4, caspase-5 and caspase-14. Cleavage of the substrate by a caspase releases pNA, which can be detected by its absorbance at 405nm.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-082-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Handling Keep cool and dry.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Caspase-1 substrate (chromogenic), Caspase-5 substrate (chromogenic)

Appearance White to yellow powder.

Formula $C_{34}H_{37}N_9O_{11}$

Formulation Lyophilized.

MW 747.7

Peptide Content 70-90%

Purity ≥98% (HPLC)

Sequence Ac-Trp-Glu-His-Asp-pNA (pNA=p-Nitroaniline)

Solubility Soluble in DMSO. Dilute with buffer, pH 7.5.

Technical Info / Product Notes λ_{max} of pNA is 400nm.

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)