

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 ₉ O 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	λmax of pNA is 400nm.

Last modified: May 29, 2024