

# Ac-WEHD-CHO

Inhibitor of caspase-1, -4, -5 and -14.

Reversible, cell permeable inhibitor of caspase-1 and -5. Also inhibits caspase-4 and -14.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-055-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Solution is stable for 1 month when stored at -20°C and protected from light.
<b>Handling</b>	Keep cool and dry.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Caspase-1 inhibitor (aldehyde), Caspase-5 inhibitor (aldehyde)
<b>Appearance</b>	White to off-white powder.
<b>Formula</b>	$C_{28}H_{33}N_7O_9$
<b>Formulation</b>	TFA salt.
<b>MW</b>	611.6
<b>Purity</b>	≥90% (HPLC)
<b>Sequence</b>	Ac-Trp-Glu-His-Asp-CHO
<b>Solubility</b>	Soluble in water (10mg/ml).

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-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)