

# Humanin (human)

## Bax inhibitor

Humanin was originally identified as an anti-apoptotic peptide of 24 aa encoded in a cDNA that rescued neuronal cells from apoptosis induced by presenilin mutants associated with familial Alzheimer's disease and by amyloid- $\beta$  protein. Mutagenesis studies showed that the cysteine at position 8 of the humanin peptide is critical for its anti-apoptotic function. It has now been shown that Bax (Bcl-2 associated X protein), an apoptosis-inducing protein, interacts with humanin. Humanin prevents the translocation of Bax from the cytosol to mitochondria. Humanin peptides also block Bax association with isolated mitochondria, and suppress cytochrome c release *in vitro*.

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-158-0110	0.5mg
ALX-158-011-M001	1mg

## 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** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** White to off-white powder.

**CAS** 330936-69-1

**Formula**  $C_{119}H_{204}N_{34}O_{32}S_2$

**MW** 2687.3

**Purity** ≥98% (HPLC)

**Sequence** H-Met-Ala-Pro-Arg-Gly-Phe-Ser-Cys-Leu-Leu-Leu-Leu-Thr-Ser-Glu-Ile-Asp-Leu-Pro-Val-Lys-Arg-Arg-Ala-OH

**Solubility** Soluble in DMSO or 5% acetic acid.

**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)