

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- β 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-158-011-MC05	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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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