

PAC-1

Caspase-3 activator

First molecule to directly activate procaspase-3 to caspase-3 *in vitro*. Has been shown to induce apoptosis in cancer cells.

Citations: 1

[View Online »](#)

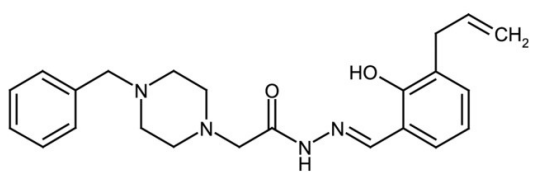
Ordering Information

[Order Online »](#)

ALX-430-146-M010	10mg
ALX-430-146-M050	50mg

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

Alternative Name 4-(Phenylmethyl)[[2-hydroxy-3-(2-propenyl)phenyl]methylene]hydrazide-1-piperazine acetic acid

Appearance White to off-white solid.

CAS 315183-21-2

Couple Target Caspase, Caspase-3

Couple Type Activator

Formula $C_{23}H_{28}N_4O_2$

MW 392.5

Purity ≥98% (HPLC (UV))

Solubility Soluble in acetonitrile, DMSO, dioxane or methanol.

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)