

Z-Asp-2,6-dichlorobenzoyloxymethylketone

Caspase inhibitor

Irreversible, cell permeable broad-spectrum caspase inhibitor.

Caspase-3 IC₅₀: 1,1µM

Caspase-6 IC₅₀: 4,1µM

Less effective on the inhibition of caspase-1- and 8-like activities.

Inhibits hypoxia-induced apoptosis in neuroblastoma cells.

Citations: 10

[View Online »](#)

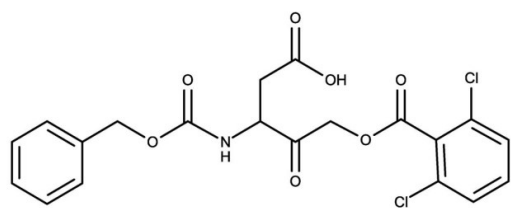
Ordering Information

[Order Online »](#)

ALX-260-029-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions should be divided into aliquots and stored at -20°C.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase family inhibitor, Z-Asp-CH ₂ -DCB, Z-L-Asp-a-[2,6-(dichlorobenzoyl)oxy]methane
Appearance	White to off-white powder.
CAS	153088-73-4
Formula	C ₂₀ H ₁₇ Cl ₂ NO ₇
MW	454.3
Purity	≥98% (TLC)
Solubility	Soluble in methanol (50mg/ml) or DMSO.

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)