

Salubrinal

Cell permeable and selective inhibitor of the phosphatase complexes that dephosphorylate eukaryotic translation initiation factor 2 subunit α (eIF-2 α). Protects ER stress-induced apoptosis ($EC_{50} \sim 15 \mu M$ in PC12 cells stimulated with 750ng/ml of tunicamycin (Prod. No. BML-CC104)).

Citations: 10

[View Online »](#)

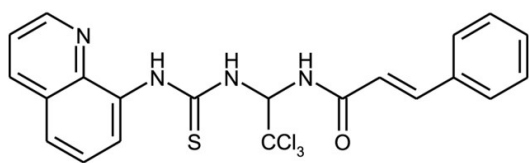
Ordering Information

[Order Online »](#)

ALX-270-428-M001	1mg
ALX-270-428-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Product may darken when exposed to air. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
CAS	405060-95-9
Formula	$C_{21}H_{17}Cl_3N_4OS$
Identity	Determined by MS, 1H-NMR and 13C-NMR.
MW	479.8
Purity	≥98% (HPLC(UV))
Solubility	Soluble in DMSO (300 mg/ml), dioxane (0.5 mg/ml), methanol (5 mg/ml) or acetonitrile (1.5 mg/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-
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)