

HA14-1

Bcl-2 ligand

Cell permeable, low molecular weight Bcl-2 ligand. Antagonizes the function of Bcl-2 and induces apoptosis. In HL-60 cells, HA4-1 (50 μ M and 4h incubation) induced apoptosis of more than 90% of the cells. Apoptosis is blocked by ZVAD-FM. IC₅₀ = ~9 μ M for competition with Flu-BakBH3 binding to Bcl-2.

Citations: 7

[View Online »](#)

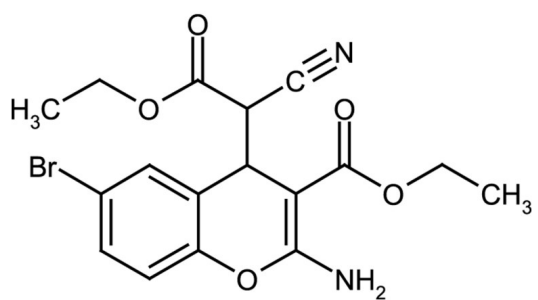
Ordering Information

[Order Online »](#)

ALX-430-101-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Unstable in solution. Prepare fresh solution for each application.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ethyl 2-amino-6-bromo-4-(1-cyano-2-ethoxy-2-oxoethyl)-4H-chromene-3-carboxylate
Appearance	Off-white to light yellow powder.
CAS	65673-63-4
Couple Target	Bcl-2
Couple Type	Ligand
Formula	$C_{17}H_{17}BrN_2O_5$
Identity	Identity determined by IR, ¹ H-NMR.
MW	409.2
Purity	≥95% (TLC)
Solubility	Soluble in DMSO (25mg/ml), dimethyl formamide or halogenated solvents; also soluble in 100% ethanol (25mg/7ml, hot).



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)