

Licochalcone A (synthetic)

Inhibitor of Bcl-2 expression

Reduces Bcl-2 protein expression.

Citations: 2

[View Online »](#)

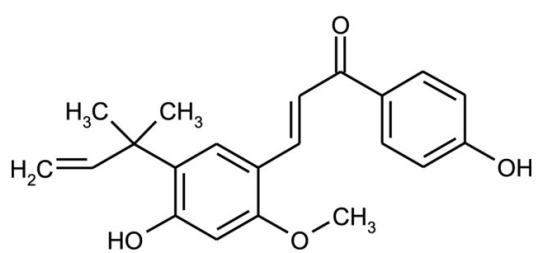
Ordering Information

[Order Online »](#)

ALX-430-124-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 6 months when stored at $\square 20^{\circ}\text{C}$. Dilute stock solution 1:10 with aqueous buffer or medium just prior to use.
Handling	Protect from light. Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C .
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(E)-3-[5-(1,1-Dimethyl-2-propenyl)-4-hydroxy-2-methoxyphenyl]-1-(4-hydroxyphenyl)-2-propen-1-one
Appearance	Yellow crystalline solid.
CAS	58749-22-7
Formula	$\text{C}_{21}\text{H}_{22}\text{O}_4$
Identity	Identity determined by NMR.
MW	338.4
Purity	$\geq 96\%$ (HPLC)
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol.

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-
be@enzolifesciences.com](mailto:info-be@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)