

Tyrphostin AG 1024

IGF-1 inhibitor

Specific inhibitor of insulin-like growth factor-1 (IGF-1) and insulin receptor kinases. Induces apoptosis by downregulation of the expression of phospho-Akt1 and Bcl-2. Also increases Bax, p53 and p21 expression.

Citations: 6

[View Online »](#)

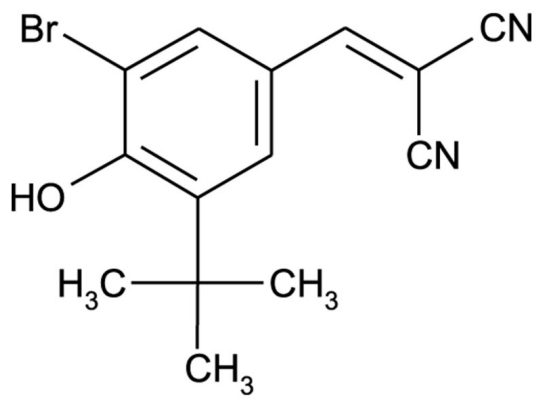
Ordering Information

[Order Online »](#)

ALX-270-217-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. In a dissolved form it should be stable for at least 1 month.
Handling	Protect from light.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Yellow powder.
CAS	65678-07-1
Couple Target	IGF
Couple Type	Inhibitor
Formula	$C_{14}H_{13}BrN_2O$
MW	305.2
Purity	≥98% (HPLC (UV))
Solubility	Soluble in acetonitrile (5mg/ml), DMSO (8mg/ml), dichloromethane (23mg/ml) or methanol (5mg/ml); insoluble in 100% ethanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com