

Damnacanthal

p56^{lck} tyrosine kinase inhibitor

Damnacanthal is a potent and selective inhibitor of p56^{lck} tyrosine kinase. It inhibits p56^{lck} autophosphorylation (IC_{50} =17 nM) as well as phosphorylation of exogenous substrates (IC_{50} =620 nM) in cell-free assays. Damnacanthal shows a 10-fold selectivity for p56^{lck} over PKA and PKC and 40-fold selectivity over four receptor tyrosine kinases as well as 7–20-fold selectivity over the homologous enzymes p60^{src} and p59^{fyn}.

Citations: 5

[View Online »](#)

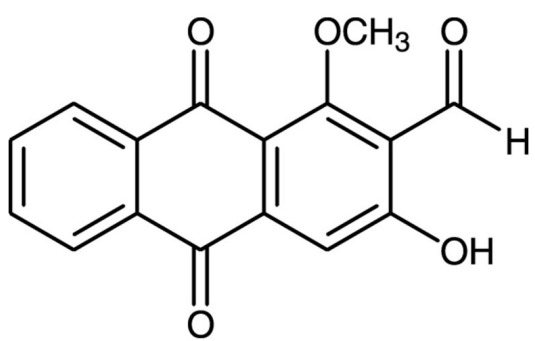
Ordering Information

[Order Online »](#)

BML-EI274-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-Hydroxy-1-methoxyanthraquinone-2-aldehyde
Appearance	Pale orange solid.
CAS	477-84-9
Couple Target	p56lck tyrosine kinase
Couple Type	Inhibitor
Formula	$C_{16}H_{10}O_5$
Identity	Determined by NMR.
MW	282.2
Purity	≥95% (HPLC)
Solubility	Soluble in 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-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)