

Daphnetin

Protein kinase inhibitor

Inhibitor of protein kinases. EGF receptor tyrosine kinase ($IC_{50}=7.87\mu M$), PKA ($IC_{50}=9.33\mu M$), PKC ($IC_{50}=25.01\mu M$). Antioxidant. Antimalarial agent.

Citations: 2

[View Online »](#)

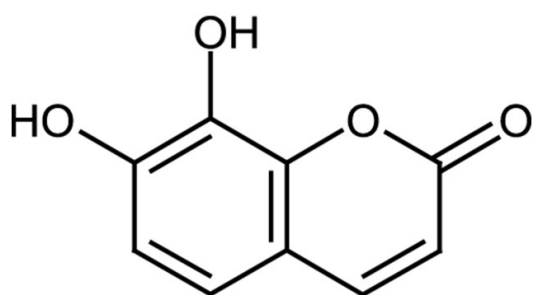
Ordering Information

[Order Online »](#)

ALX-270-281-M005	5mg
ALX-270-281-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 7,8-Dihydroxycoumarin

Appearance White to rose powder.

CAS 486-35-1

Couple Target EGFR, PKA, PKC

Couple Type Inhibitor

Formula $C_9H_6O_4$

Identity Determined by NMR.

MW 178.1

Purity $\geq 97\%$ (HPLC)

RTECS GN6382600

Solubility Soluble in 100% ethanol, methanol or dimethyl formamide.

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)