

Neobavaisoflavone

DNA polymerase inhibitor

Inhibits platelet aggregation. DNA polymerase inhibitor. Shows antifungal activity.

Citations: 5

[View Online »](#)

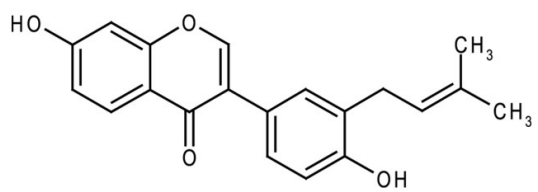
Ordering Information

[Order Online »](#)

ALX-350-146-M001	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

Appearance	Yellow solid.
CAS	41060-15-5
Couple Target	DNA polymerase
Couple Type	Inhibitor
Formula	$C_{20}H_{18}O_4$
Identity	Determined by 1H-NMR and MS.
MW	322.4
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol.
Source	Isolated from plant <i>Psoralea corylifolia</i> .

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)