

Genistein (synthetic)

Tyrosine kinase inhibitor

Tyrosine protein kinase inhibitor. Inhibits phosphorylation of EGFR kinase. Inhibits tumor cell proliferation and induces tumor cell differentiation. Inhibits topoisomerase II activity *in vivo*. Produces cell cycle arrest and apoptosis. Potent inhibitor of α -glucosidase, $IC_{50} = 50\text{nM}$. Direct inhibitor of insulin-induced glucose uptake in adipocytes ($IC_{50} = 20\mu\text{M}$).

Citations: 26

[View Online »](#)

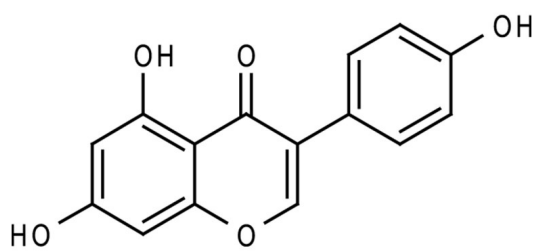
Ordering Information

[Order Online »](#)

ALX-350-006-M010	10mg
ALX-350-006-M025	25mg
ALX-350-006-M050	50mg

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	4',5,7-Trihydroxyisoflavone
Appearance	Creamy off-white solid.
CAS	446-72-0
Couple Target	Glucosidase, Topoisomerase
Couple Type	Inhibitor
Formula	$C_{15}H_{10}O_5$
MI	14: 4391
MW	270.2
Purity	≥98% (HPLC)
RTECS	NR2392000
Solubility	Soluble in DMSO (30mg/ml), 100% ethanol (4mg/ml) or dilute aqueous base; insoluble in water.
Source	Synthetic.

Technical Info / Product Notes



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

Replacement for ADI-HPK-103

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)

Last modified: May 29, 2024