

Rottlerin

PKCδ inhibitor

Mitochondrial uncoupler that depolarizes the mitochondrial membrane potential, reduces cellular ATP levels, activates 5'-AMP-activated protein kinase (AMPK) and affects mitochondrial production of reactive oxygen species (ROS). Potent activator of multiple Ca^{2+} -sensitive K^{+} channels. Blocks several kinases and non-kinase proteins *in vitro*. Has been widely-used as a selective inhibitor of protein kinase Cδ (PKCδ) (IC_{50} =3-6μM). Although there is extensive published documentation to support the use of rottlerin as a selective PKCδ inhibitor, there has been some controversy in the literature over this claim. Some of this controversy may arise from the fact that rottlerin is a promiscuous inhibitor and therefore may not always display the same properties as would be expected from a classical 1:1 inhibitor. However, there are studies indicate that rottlerin has no direct effect on PKCδ and that it should not be used to determine the involvement of PKCδ in biological processes. Induces autophagy by inhibition of mTORC1 signaling.

Citations: 11

[View Online »](#)

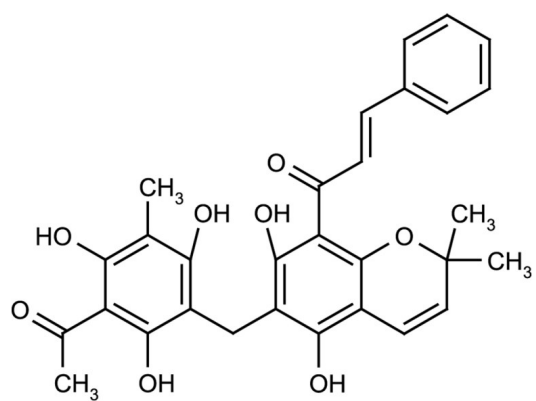
Ordering Information

[Order Online »](#)

ALX-350-075-M010	10mg
ALX-350-075-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 Mallotoxin

Appearance Orange to brown solid.

CAS 82-08-6

Couple Target PKC

Couple Type Inhibitor

Formula $C_{30}H_{28}O_8$

MI 14: 8272

MW 516.5

Purity ≥98% (HPLC)

RTECS AM6913800

Solubility Soluble in DMSO (up to 50mg/ml).

Source Isolated from *Mallotus philippinensis*.



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)

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)