

Rottlerin

PKC δ 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 Ca²⁺-sensitive K⁺ channels. Blocks several kinases and non-kinase proteins *in vitro*. Has been widely-used as a selective inhibitor of protein kinase C δ (PKC δ) (IC₅₀=3-6 μ M). Although there is extensive published documentation to support the use of rottlerin as a selective PKC δ 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 PKC δ and that it should not be used to determine the involvement of PKC δ in biological processes. Induces autophagy by inhibition of mTORC1 signaling.

Citations: 11

[View Online »](#)

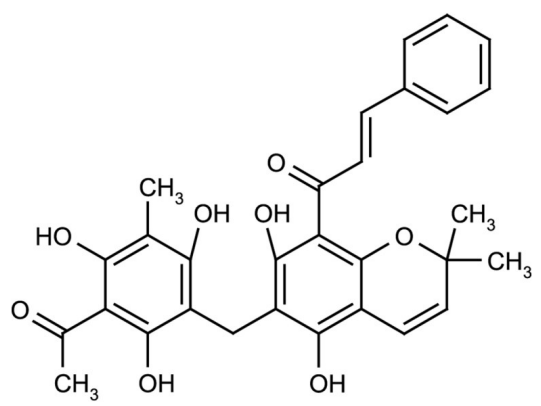
Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-350-075-M010 | 10mg |
|------------------|------|

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 <i>Mallotus philippinensis</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-
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)