

Pterostilbene

CYP1A1 inhibitor

Cell permeable natural methoxylated analog of resveratrol (Prod. No. BML-FR104). Antioxidant, antiproliferative, anti-inflammatory, anti-hyperglycemic and anti-diabetic agent. Induces apoptosis. Induces Autophagy. Very potent inhibitor of cytochrome P450 1A1 (CYP1A1) with $K_i=0.57\mu\text{M}$. Inhibits activator protein 1 (AP-1) and NF- κ B activation. Moderately inhibits COX-1 and COX-2 ($\text{IC}_{50}=19.8\mu\text{M}$ and $83.9\mu\text{M}$).

Citations: 22

[View Online »](#)

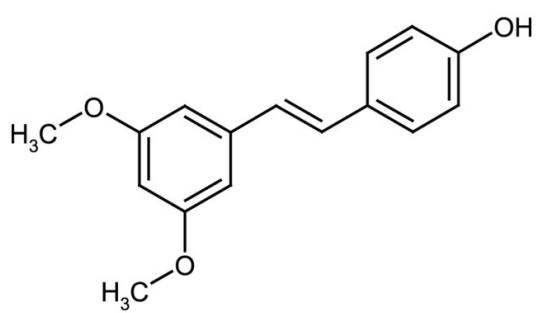
Ordering Information

[Order Online »](#)

ALX-385-034-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and oxygen.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	trans-3,5-Dimethoxy-4'-hydroxystilbene
Appearance	Off-white to light brown powder.
CAS	537-42-8
Couple Target	COX, Jun
Couple Type	Inhibitor
Formula	$C_{16}H_{16}O_3$
Identity	Identity determined by MS and 1H -NMR.
MW	256.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (20mg/ml) or methanol; insoluble in water.
Source	Synthetic.

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)