

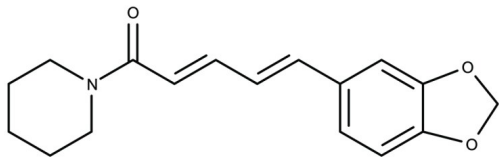
Piperine

CYP3A4 inhibitor. Natural product.

Piperine is a natural product, isolated from black pepper, with many biological effects. It increases the metabolic rate of fast twitch skeletal fibers, via destabilization of the super-relaxed state of myosin heads, without effecting slow twitch or cardiac fibers. TRPV1 agonist (EC50 = 38 μ M), and inhibits human P-glycoprotein CYP3A4 and UDP-glucose dehydrogenase and glucuronidation.

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored unopened at -20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white solid.
CAS	94-62-2
Couple Target	Cytochrome P450, TRP channel
Couple Type	Activator, Inhibitor
Formula	$C_{17}H_{19}NO_3$
Formulation	Lyophilized.
Identity	Identity determined by NMR.
MW	285.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25 mg/ml).

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)