

# 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 (EC<sub>50</sub> = 38  $\mu$ M), and inhibits human P-glycoprotein CYP3A4 and UDP-glucose dehydrogenase and glucuronidation.

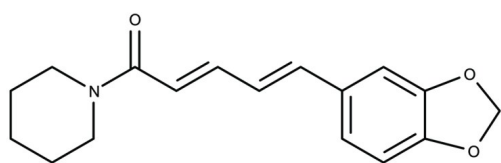
## Ordering Information

[Order Online »](#)

ENZ-CHM268-0250	250mg
-----------------	-------

## 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).



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)