

# Resveratrol

**Antioxidant. SIRT1 activator.**

Phenolic antioxidant found in grapes and wine. Potent activator of human deacetylase SIRT1 (maximum effect at 100-200 $\mu$ M). Increases longevity in yeast. Displays antiproliferative effect on human breast epithelial cells. Displays a variety of pharmacological activities including: estrogen receptor agonist activity, inhibition of ribonucleotide reductase and DNA synthesis, inhibition of eicosanoid synthesis, cell cycle arrest at S/G2 phase and induction of apoptosis. Inhibits PKD *in vitro* (IC<sub>50</sub>=200 $\mu$ M) and *in vivo* (IC<sub>50</sub>=800 $\mu$ M), but not PKC isoforms. Protects against 4-hydroxynonenal (4-HNE) induced oxidative stress and apoptosis. Shows cancer chemopreventive activity. Specific inhibitor of cyclooxygenase-1 (COX-1). Inhibits the hydroperoxidase activity of COX-1.

Citations: 38

[View Online »](#)

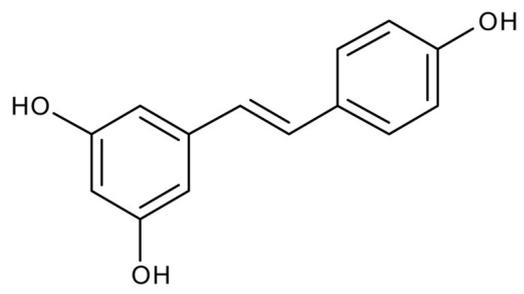
## Ordering Information

[Order Online »](#)

BML-FR104-0100	100mg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	trans-3,4',5-Trihydroxystilbene
<b>Appearance</b>	White to yellow solid.
<b>CAS</b>	501-36-0
<b>Couple Target</b>	COX, Estrogen receptor, PKD, SIRT, SIRT1
<b>Couple Type</b>	Activator, Inhibitor, Ligand
<b>Formula</b>	$C_{14}H_{12}O_3$
<b>MI</b>	<b>14:</b> 8158
<b>MW</b>	228.2
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in DMSO (65mg/ml), 100% ethanol or dimethyl formamide; slightly soluble in water (0.1mg/ml).
<b>Technical Info / Product Notes</b>	Note: Product is not sterile.

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)