

Resveratrol

Antioxidant. SIRT1 activator.

Phenolic antioxidant found in grapes and wine. Potent activator of human deacetylase SIRT1 (maximum effect at 100-200 μ 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* (IC₅₀=200 μ M) and *in vivo* (IC₅₀=800 μ 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 »](#)

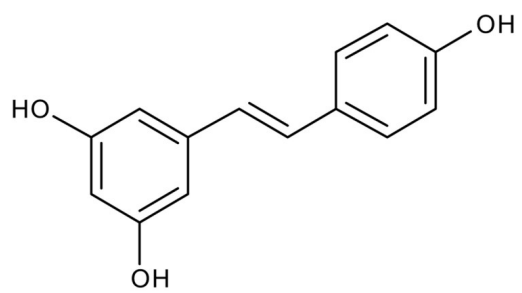
Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	trans-3,4',5-Trihydroxystilbene
Appearance	White to yellow solid.
CAS	501-36-0
Couple Target	COX, Estrogen receptor, PKD, SIRT, SIRT1
Couple Type	Activator, Inhibitor, Ligand
Formula	$C_{14}H_{12}O_3$
MI	14: 8158
MW	228.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (65mg/ml), 100% ethanol or dimethyl formamide; slightly soluble in water (0.1mg/ml).

Technical Info / Product Notes Note: Product is not sterile.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com