

Butein

Tyrosine kinase inhibitor

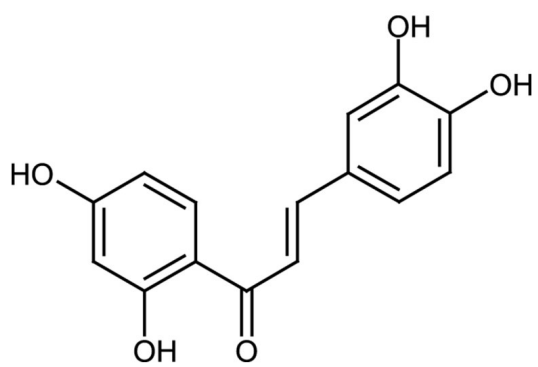
Plant polyphenol. Specific tyrosine kinase inhibitor. Potently inhibits the tyrosine kinase activity of the EGF receptor and p60^{c-src}. Potent antioxidant and anti-inflammatory agent. Inhibits glutathione reductase and rat liver glutathione S-transferase. Activator of human deacetylase SIRT1. Inhibits aromatase, showing chemopreventive properties. Directly inhibits IKK.

Citations: 15

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep cool and dry under inert gas. Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2',3,4,4'-Tetrahydroxychalcone
Appearance	Yellow solid.
CAS	487-52-5
Couple Target	Cytochrome P450, EGFR, Glutathione reductase, GST, IKK, SIRT, SIRT1
Couple Type	Activator, Inhibitor
Formula	$C_{15}H_{12}O_5$
Identity	Identity determined by 1H-NMR, 13C-NMR and MS.
MW	272.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>50mg/ml) or methanol; insoluble in water.

Last modified: May 29, 2024



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