

Cinnamtannin B-1

Potent antioxidant

A-type proanthocyanidin contained in several plant species such as *Laurus nobilis* L. , *Vaccinium vitis-idaea*, *Parameria laevigata*, *Cinnamomum zeylanicum*, *Lindera umbellata*, *Metaxya rostrata*. Potent antioxidant, TEAC method (nM Trolox) = 0.95. Protective agent against oxidative stress and apoptosis in human platelets.

Inhibitory activity against cyclooxygenase-2 (COX-2) *in vitro*.

DPPH radical scavenging IC₅₀ 6.14μM

Lipid peroxidation inhibition 159.74

Nitric oxide inhibition 10.28

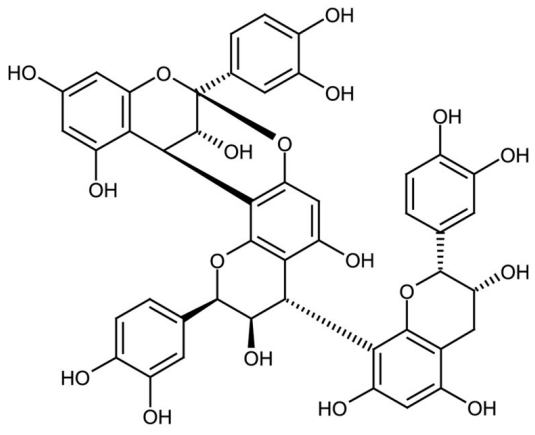
Antibacterial against *B.subtilis* (9mm, 20mg/ml)

Citations: 18

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Epicatechin-(4β→8,2β→O→7)-epicatechin-(4α→8)-epicatechin
Appearance	Brown solid.
CAS	88082-60-4
Couple Target	COX
Couple Type	Inhibitor
Formula	$C_{45}H_{36}O_{18}$
Identity	Determined by 1H-NMR.
MW	864.7
Purity	≥95% (HPLC)
Solubility	Soluble in 100% ethanol, methanol or in a mixture of DMSO and water.
Source	Isolated from <i>Laurus nobilis</i> L.

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