

# (-)-Epigallocatechin gallate

Inhibitor of NOS, telomerase and Dnmt

Antitumor reagent. Antioxidant. Protects cells from lipid peroxidation and DNA damage induced by reactive free radicals. Inhibits inducible nitric oxide synthase (iNOS; NOS II). Chemopreventive anticancer agent. Induces apoptosis in human cancer cell lines. Inhibits MAP kinase mediated signalling pathways. Inhibits angiogenesis. Inhibits telomerase and DNA methyltransferase. Anti-inflammatory agent.

Citations: 36

[View Online »](#)

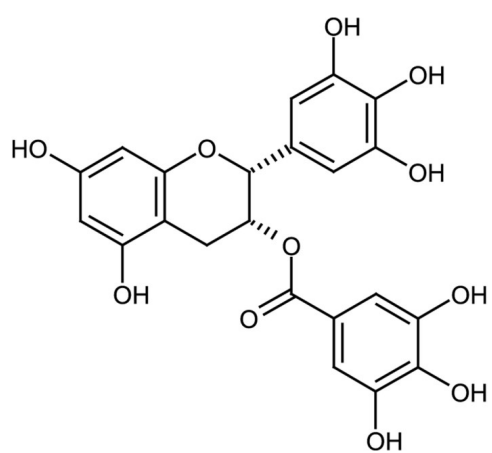
## Ordering Information

[Order Online »](#)

ALX-270-263-M010	10mg
ALX-270-263-M050	50mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO are stable for up to 3 months when stored at -20°C.
<b>Handling</b>	Protect from light. Keep under inert gas. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	EGCG, (2R,3R)-2-(3,4,5-Trihydroxyphenyl)-3,4-dihydro-1[2H]-benzopyran-3,5,7-triol 3-(3,4,5-trihydroxybenzoate)
<b>Appearance</b>	White powder.
<b>CAS</b>	989-51-5
<b>Couple Target</b>	Dnmt, NOS, Telomerase
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{22}H_{18}O_{11}$
<b>MI</b>	14: 3526
<b>MW</b>	458.4
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in 100% ethanol, dimethyl formamide or DMSO; slightly soluble in water.
<b>Source</b>	Isolated from green tea.



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)