

Gallotannin

Inhibitor of PARG and eNOS

Inhibitor of poly(ADP-ribose) glycohydrolase (PARG). Cytoprotective in oxidatively stressed cells. Inhibitor of endothelial nitric oxide synthase (eNOS; NOS III) and weak inhibitor of inducible (iNOS; NOS II) and neuronal nitric oxide synthase (nNOS; NOS I). Induces cyclooxygenase-2 (COX-2) expression. Free radical scavenger.

Citations: 12

[View Online »](#)

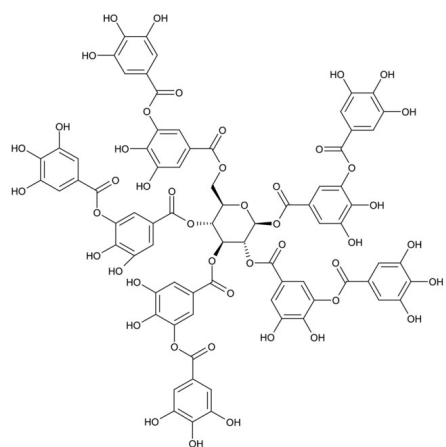
Ordering Information

[Order Online »](#)

ALX-270-418-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tannic acid

Appearance Beige to light brown powder.

CAS 1401-55-4

Couple Target NOS, PARG

Couple Type Inhibitor

Formula $C_{76}H_{52}O_{46}$

MI 14: 9052

MW 1701.2

Purity ≥95%

RTECS WW5075000

Solubility Soluble in water, 100% ethanol, acetone or warm glycerol.

Source Occurs in the bark and fruit of many plants, notably in the bark of the oak species, in sumac and myrobalan.



ENZO LIFE SCIENCES,
INC.
Phone: +1 909 440 655
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