

Koningic acid

Glyceraldehyde-3-phosphate dehydrogenase inhibitor

Koningic acid (KA) is a potent and selective inhibitor of glyceraldehyde-3-phosphate dehydrogenase (GAPDH). Inhibition is irreversible and proceeds via nucleophilic attack of an active site cysteine on the epoxide moiety, for rabbit muscle GAPDH3. KA can selectively kill high-glycolytic cancer cells via glucose dependent ATP depletion. Has been used in a predictive model for selective targeting of the Warburg effect, the most prominent hallmark of cancer cell metabolism.

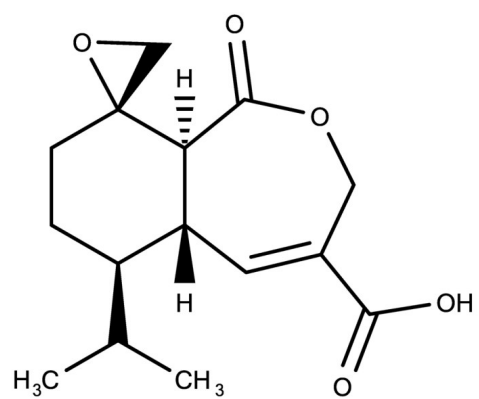
Ordering Information

[Order Online »](#)

ENZ-CHM494-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (2'S,5aS,6R,9aS)-1,5a,6,7,8,9a-Hexahydro-6-(1-methylethyl)-1-oxo-spiro[2-benzoxepin-9(3H)2'-oxirane]-4-carboxylic acid

Appearance Off-white solid.

CAS 74310-84-2

Couple Target GAPDH

Couple Type Inhibitor

Formula $C_{15}H_{20}O_5$

Identity Determined by NMR.

MW 280.32

Purity ≥95% (HPLC)

Solubility Soluble in water (1 mg/ml).



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-fr@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