

# 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.

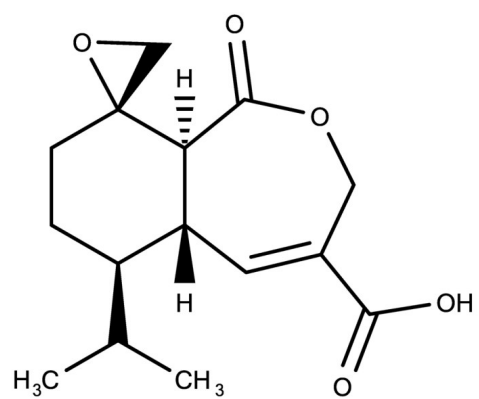
### Ordering Information

[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<sub>15</sub>H<sub>20</sub>O<sub>5</sub>

**Identity** Determined by NMR.

**MW** 280.32

**Purity** ≥95% (HPLC)

**Solubility** Soluble in water (1 mg/ml).

**Last modified: February 27, 2025**



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)