

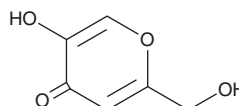
PRODUCT INFORMATION



Kojic Acid

Item No. 22712

CAS Registry No.: 501-30-4
Formal Name: 5-hydroxy-2-(hydroxymethyl)-4H-pyran-4-one
Synonym: NSC 1942
MF: $C_6H_6O_4$
FW: 142.1
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 218, 272 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Kojic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the kojic acid in the solvent of choice. Kojic acid is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. The solubility of kojic acid in these solvents is approximately 5 and 16 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of kojic acid can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of kojic acid in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Kojic acid is a fungal metabolite that inhibits tyrosinase, an enzyme involved in melanin synthesis, with an IC_{50} value of $30.6 \mu\text{M}$ for mushroom tyrosinase.¹ It decreases growth of *Leishmania* parasites *in vitro* and *in vivo*, and it protects mice from damage induced by gamma irradiation.^{2,3} Kojic acid has been used as a food additive to prevent enzymatic browning.⁴ Formulations containing kojic acid are used in the treatment of hyperpigmentation disorders.⁵

References

1. Lee, Y.S., Park, J.H., Kim, M.H., *et al.* Synthesis of tyrosinase inhibitory kojic acid derivative. *Arch. Pharm. (Weinheim)* **339**(3), 111-114 (2006).
2. Rodrigues, A.P., Farias, L.H., Carvalho, A.S., *et al.* A novel function for kojic acid, a secondary metabolite from *Aspergillus* fungi, as antileishmanial agent. *PLoS One* **9**(3), e91259 (2014).
3. Wang, K., Liu, C., Di, C.-J., *et al.* Kojic acid protects C57BL/6 mice from gamma-irradiation induced damage. *Asian Pac. J. Cancer Prev.* **15**(1), 291-297 (2014).
4. Burdock, G.A., Soni, M.G., and Carabin, I.G. Evaluation of health aspects of kojic acid in food. *Regul. Toxicol. Pharmacol.* **33**(1), 80-101 (2001).
5. Kim, Y.M., Yun, J., Lee, C.K., *et al.* Oxyresveratrol and hydroxystilbene compounds. Inhibitory effect on tyrosinase and mechanism of action. *J. Biol. Chem.* **277**(18), 16340-16344 (2002).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/22/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM