

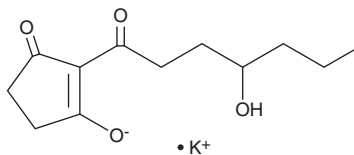
PRODUCT INFORMATION



CAY10799 (potassium salt)

Item No. 40227

CAS Registry No.: 34200-97-0
Formal Name: 3-hydroxy-2-(4-hydroxy-1-oxoheptyl)-2-cyclopenten-1-one, monopotassium salt
MF: C₁₂H₁₇O₄ • K
FW: 264.4
Purity: ≥98%
Supplied as: A powder
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

CAY10799 (potassium salt) is supplied as a powder. A stock solution may be made by dissolving the CAY10799 (potassium salt) in the solvent of choice, which should be purged with an inert gas. CAY10799 (potassium salt) is soluble in the organic solvent DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10799 is a derivative of the fungal metabolite and tyrosine hydroxylase inhibitor oudenone.^{1,2} It undergoes pH-dependent spontaneous cyclization to oudenone in aqueous solutions.¹

References

1. Tsantrizos, Y.S., Zhou, F., Famili, P., *et al.* Biosynthesis of the hypotensive metabolite oudenone by *Oudemansiella radicata*. 1. Intact incorporation of a tetraketide chain elongation intermediate. *J. Org. Chem.* **60**(21), 6922-6929 (1995).
2. Koizumi, S., Nagatsu, T., Iinuma, H., *et al.* Inhibition of phenylalanine hydroxylase, a pterin-requiring monooxygenase, by oudenone and its derivatives. *J. Antibiot. (Tokyo)* **35**(4), 458-462 (1982).

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, 11/14/2023

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