

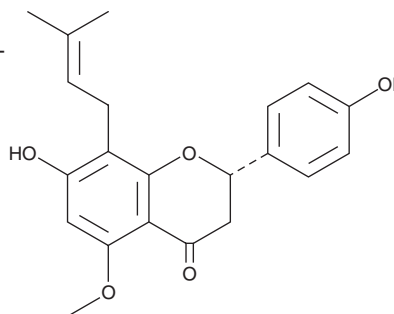
PRODUCT INFORMATION



(-)-Isoxanthohumol

Item No. 25105

CAS Registry No.: 70872-29-6
Formal Name: (2S)-2,3-dihydro-7-hydroxy-2-(4-hydroxyphenyl)-5-methoxy-8-(3-methyl-2-buten-1-yl)-4H-1-benzopyran-4-one
MF: C₂₁H₂₂O₅
FW: 354.4
Purity: ≥98%
UV/Vis.: λ_{max}: 290 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

(-)-Isoxanthohumol is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-isoxanthohumol in the solvent of choice. (-)-Isoxanthohumol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of (-)-isoxanthohumol in ethanol and DMF is approximately 3 mg/ml and approximately 2 mg/ml in DMSO.

(-)-Isoxanthohumol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (-)-isoxanthohumol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (-)-Isoxanthohumol has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(-)-Isoxanthohumol is an enantiomer of isoxanthohumol (Item No. 23233). It is found in similar amounts as (+)-isoxanthohumol in beer, where xanthohumol is converted to isoxanthohumol during the brewing process.¹

Reference

1. Kodama, S., Yamamoto, A., Sato, A., *et al.* Enantioseparation of isoxanthohumol in beer by hydroxypropyl-gamma-cyclodextrin-modified micellar electrokinetic chromatography. *J. Agric. Food Chem.* **55**(16), 6547-6552 (2007).

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, 12/20/2022

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