

PRODUCT INFORMATION



Hesperidin methylchalcone/*C. aurantium* L. Extract (hydrate)

Item No. 30884

Formal Name: (2E)-1-[4-[[6-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranosyl]oxy]-2-hydroxy-6-methoxyphenyl]-3-(3-hydroxy-4-methoxyphenyl)-2-propen-1-one, hydrate

MF: $C_{29}H_{36}O_{15} \cdot XH_2O$

FW: 624.6

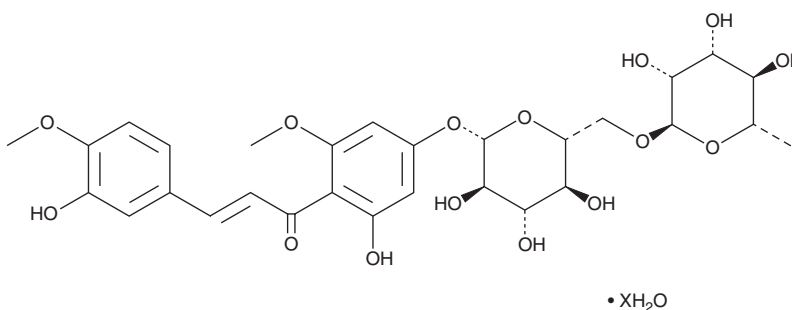
UV/Vis.: λ_{max} : 284 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

Item Origin: Plant/*Citrus aurantium* L.



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

Laboratory Procedures

Hesperidin methylchalcone/*C. aurantium* L. extract (hydrate) is supplied as a solid. A stock solution may be made by dissolving the hesperidin methylchalcone/*C. aurantium* L. extract (hydrate) in the solvent of choice, which should be purged with an inert gas. Hesperidin methylchalcone/*C. aurantium* L. extract (hydrate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of hesperidin methylchalcone/*C. aurantium* L. extract (hydrate) in these solvents is approximately 20 mg/ml.

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 hesperidin methylchalcone/*C. aurantium* L. extract (hydrate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of hesperidin methylchalcone/*C. aurantium* L. extract (hydrate) in PBS, pH 7.2, is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Hesperidin methylchalcone/*C. aurantium* L. extract is a *C. aurantium* extract that contains hesperidin methylchalcone, among other compounds.

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, 11/08/2022

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