

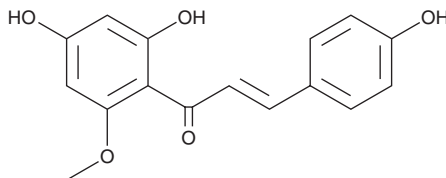
PRODUCT INFORMATION



Helichrysetin

Item No. 43168

CAS Registry No.: 62014-87-3
Formal Name: (2E)-1-(2,4-dihydroxy-6-methoxyphenyl)-3-(4-hydroxyphenyl)-2-propen-1-one
Synonym: 6'-Methoxy-2',4,4'-trihydroxy-chalcone
MF: $C_{16}H_{14}O_5$
FW: 286.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Xerochrysum bracteatum*



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

Laboratory Procedures

Helichrysetin is supplied as a solid. A stock solution may be made by dissolving the helichrysetin in the solvent of choice, which should be purged with an inert gas. Helichrysetin is soluble (≥ 10 mg/ml) in ethanol.

Description

Helichrysetin is a chalcone that has been found in *H. cymosum* and has diverse biological activities.¹⁻⁴ It inhibits α -glucosidase and reduces ferric to ferrous iron in a ferric reducing antioxidant power (FRAP) assay (IC_{50} s = 18.16 and 279.93 μM , respectively).¹ Helichrysetin (1 mg/ml) is active against *B. subtilis*.² It also inhibits Middle East respiratory syndrome coronavirus (MERS-CoV) main protease (M^{pro}), also known as 3C-like protease ($3CL^{pro}$; IC_{50} = 67.04 μM).³ *In vivo*, helichrysetin (10 mg/kg) reduces tumor growth in an MCF10DCIS.com breast cancer mouse xenograft model.⁴

References

1. Jadalla, B.M.I.S., Moser, J.J., Sharma, R., *et al.* In vitro alpha-glucosidase and alpha-amylase inhibitory activities and antioxidant capacity of *Helichrysum cymosum* and *Helichrysum pandurifolium* schrank constituents. *Separations* **9**(8), 190 (2022).
2. Malolo, F.A., Bissoue Noug, A., Kakam, A., *et al.* Protease-inhibiting, molecular modeling and antimicrobial activities of extracts and constituents from *Helichrysum foetidum* and *Helichrysum mechowianum* (compositae). *Chem. Cent. J.* **9**, 32 (2015).
3. Jo, S., Kim, H., Kim, S., *et al.* Characteristics of flavonoids as potent MERS-CoV 3C-like protease inhibitors. *Chem. Biol. Drug Des.* **94**(6), 2023-2030 (2019).
4. Liu, Y., Pandey, P.R., Sharma, S., *et al.* ID2 and GJB2 promote early-stage breast cancer progression by regulating cancer stemness. *Breast Cancer Res. Treat.* **175**(1), 77-90 (2019).

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, 04/16/2025

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