

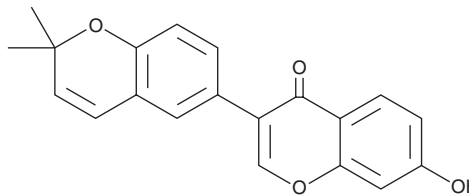
PRODUCT INFORMATION



Corylin

Item No. 45980

CAS Registry No.: 53947-92-5
Formal Name: 3-(2,2-dimethyl-2H-1-benzopyran-6-yl)-7-hydroxy-4H-1-benzopyran-4-one
MF: $C_{20}H_{16}O_4$
FW: 320.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: 20°C
Stability: ≥ 4 years
Item Origin: Plant/*Psoralea corylifolia*



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

Laboratory Procedures

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

Description

Corylin is a flavonoid that has been found in *P. corylifolia* and has diverse biological activities.¹⁻⁸ It inhibits carboxylesterase 1 (CES1; $K_i = 0.7 \mu\text{M}$), aldose reductase ($\text{IC}_{50} = 7.5 \mu\text{M}$), and IL-6-induced STAT3 activation in a reporter assay using Hep3B cells ($\text{IC}_{50} = 1.37 \mu\text{M}$).¹⁻³ Corylin ($10 \mu\text{M}$) inhibits the migration and invasion of MCF-7 and MDA-MB-231 breast cancer cells and LPS-induced prostaglandin E_2 (PGE_2 ; Item No. 14010) and nitric oxide (NO) production in RAW 264.7 cells.^{4,5} It induces osteogenesis in micromass cultures from rat E13 limb bud cells.⁶ Corylin decreases body weight and white adipose tissue weight and induces adipocyte browning in a mouse model of diet-induced obesity and increases survival in aged mice fed a high-fat diet.^{7,8}

References

1. Sun, D.X., Ge, G.B., Dong, P.P., et al. Inhibition behavior of fructus psoraleae's ingredients towards human carboxylesterase 1 (hCES1). *Xenobiotica* **46**(6), 503-510 (2016).
2. Shim, S.H., Kim, Y., Lee, J.Y., et al. Aldose reductase inhibitory activity of the compounds from the seed of *Psoralea corylifolia*. *J. Korean Soc. Appl. Biol. Chem.* **52**(5), 568-572 (2009).
3. Lee, S.W., Yun, B.R., Kim, M.H., et al. Phenolic compounds isolated from *Psoralea corylifolia* inhibit IL-6-induced STAT3 activation. *Planta Med.* **78**(9), 903-906 (2012).
4. Liu, S., Wang, L., and Zhang, R. Corylin suppresses metastasis of breast cancer cells by modulating miR-34c/LINC00963 target. *Libyan J. Med.* **16**(1), 1883224 (2021).
5. Hung, Y.L., Fang, S.H., Wang, S.C., et al. Corylin protects LPS-induced sepsis and attenuates LPS-induced inflammatory response. *Sci. Rep.* **7**, 46299 (2017).
6. Yu, A.X., Xu, M.L., Yao, P., et al. Corylin, a flavonoid derived from *Psoralea Fructus*, induces osteoblastic differentiation via estrogen and Wnt/ β -catenin signaling pathways. *FASEB J.* **34**(3), 4311-4328 (2020).
7. Chen, C.C., Kuo, C.H., Leu, Y.L., et al. Corylin reduces obesity and insulin resistance and promotes adipose tissue browning through SIRT-1 and $\beta 3$ -AR activation. *Pharmacol. Res.* **164**, 105291 (78370).
8. Wang, T.H., Tseng, W.C., Leu, Y.L., et al. The flavonoid corylin exhibits lifespan extension properties in mouse. *Nat. Commun.* **13**(1), 1238 (2022).

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, 07/15/2026

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