

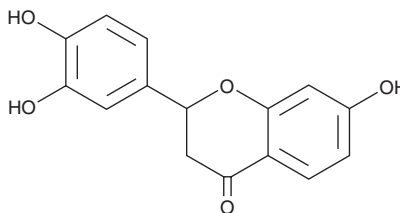
PRODUCT INFORMATION



Butin

Item No. 40268

CAS Registry No.: 21913-99-5
Formal Name: 2-(3,4-dihydroxyphenyl)-2,3-dihydro-7-hydroxy-4H-1-benzopyran-4-one
Synonym: 3',4',7-Trihydroxyflavanone
MF: $C_{15}H_{12}O_5$
FW: 272.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Butin is supplied as a solid. A stock solution may be made by dissolving the butin in the solvent of choice, which should be purged with an inert gas. Butin is soluble in acetonitrile and methanol.

Description

Butin is a flavanone that has been found in *B. monosperma* and has diverse biological activities.¹⁻³ It induces melanogenesis in B16/F10 melanoma cells when used at a concentration of $1\text{ }\mu\text{M}$.¹ Butin ($1\text{--}4\text{ }\mu\text{g/ml}$) decreases DNA damage induced by the actin polymerization inhibitor cytochalasin B (Item No. 11328) in isolated human peripheral blood lymphocytes in a cytokinesis-block micronucleus (CBMN) assay.² *In vivo*, butin ($5\text{--}20\text{ mg/animal}$) inhibits implantation and decreases the number of pups when administered to pregnant rats.³ It increases uterine weight in ovariectomized rats when administered at a dose of 0.5 mg/animal .

References

1. Lv, J., Zha, X., Pang, S., et al. Synthesis and melanogenesis evaluation of 3',4',7-trihydroxyflavanone derivatives and characterization of flavanone-BODIPY. *Bioorg. Med. Chem. Lett.* **25(7)**, 1607-1610 (2015).
2. Milosavljevic, S., Djordjevic, I., Mandic, B., et al. Flavonoids of the heartwood of *Cotinus coggygria* Scop. showing protective effect on human lymphocyte DNA. *Nat. Prod. Commun.* **16(12)**, 1-8 (2021).
3. Bhargava, S.K. Estrogenic and postcoital contraceptive activity in rats of butin isolated from *Butea monosperma* seed. *J. Ethnopharmacol.* **18(1)**, 95-101 (1986).

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, 04/23/2024

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