

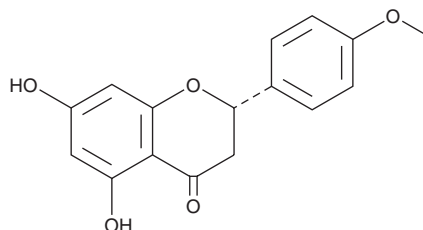
PRODUCT INFORMATION



Isosakuranetin

Item No. 25198

CAS Registry No.: 480-43-3
Formal Name: (2S)-2,3-dihydro-5,7-dihydroxy-2-(4-methoxyphenyl)-4H-1-benzopyran-4-one
MF: C₁₆H₁₄O₅
FW: 286.3
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 226, 289 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

Isosakuranetin is supplied as a crystalline solid. A stock solution may be made by dissolving the isosakuranetin in the solvent of choice, which should be purged with an inert gas. Isosakuranetin is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of isosakuranetin in these solvents is approximately 25 and 30 mg/ml, respectively.

Isosakuranetin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, isosakuranetin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Isosakuranetin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Isosakuranetin is a flavanone that has been found in *Citrus* species and has diverse biological activities.¹⁻⁷ It inhibits calcium uptake induced by pregnenolone sulfate (PregS; Item No. 21004) in HEK293 cells expressing mouse transient receptor potential melastatin 3 (TRPM3; IC₅₀ = 50 nM).¹ Isosakuranetin is selective for TRPM3 over TRPM1, TRPM8, and TRP vanilloid-related 1 (TRPV1) when used at a concentration of 10 μM. *In vivo*, isosakuranetin (2 mg/kg) increases latency to first pain-related response in mice in a hot plate test and reduces the number and duration of PregS-induced nocifensive responses, such as paw licking, shaking, or lifting, in mice. It also reduces systolic blood pressure in spontaneously hypertensive rats when administered at a dose of 10 mg/kg.² Isosakuranetin (20 μM) inhibits UV-B-induced matrix metalloproteinase-1 (MMP-1) expression by 90% in HaCaT human keratinocyte cells, as well as phosphorylation of ERK1/2 when used at a concentration of 50 μM.³ It blocks hydrogen peroxide-induced increases in reactive oxygen species (ROS), intracellular calcium concentration, caspase-3 activity, and JNK phosphorylation, as well as decreases in catalase activity, in PC12 cells when used at a concentration of 0.8 μM.⁴ Isosakuranetin also has antibacterial and trypanocidal activity against *M. tuberculosis*, *C. neoformans*, and *T. cruzi*.⁵⁻⁷

References

1. Straub, I., Krügel, U., Mohr, F., et al. *Mol. Pharmacol.* **84**(5), 736-750 (2013).
2. Maruyama, H., Sumitou, Y., Sakamoto, T., et al. *Biol. Pharm. Bull.* **32**(7), 1244-1250 (2009).
3. Jung, H., Lee, E.H., Lee, T.H., et al. *Int. J. Mol. Sci.* **17**(9), E1449 (2016).
4. Hwang, S.-L. and Yen, G.-C. *J. Agric. Food Chem.* **57**(6), 2576-2582 (2009).
5. Suksamrarn, A., Chotipong, A., Suavansri, T., et al. *Arch. Pharm. Res.* **27**(5), 507-511
6. da Silva Filho, A.A., de Sousa, J.P., Soares, S., et al. *Z. Naturforsch. C.* **63**(1-2), 40-46 (2008).
7. da Silva Filho, A.A., Pires Bueno, P.C., Gregório, L.E., et al. *J. Pharm. Pharmacol.* **56**(9), 1195-1199 (2004).

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, 12/02/2022

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