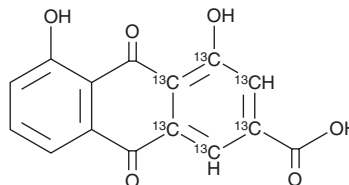


PRODUCT INFORMATION



Rhein-¹³C₆
Item No. 42286

Formal Name: 9,10-dihydro-4,5-dihydroxy-9,10-dioxo-2-anthracenecarboxylic acid-1,2,3,4,4a,9a-¹³C₆
Synonym: Rheic Acid-¹³C₆
MF: C₉[¹³C]₆H₈O₆
FW: 290.2
Purity: ≥95%
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

Rhein-¹³C₆ is supplied as a solid. A stock solution may be made by dissolving the rhein-¹³C₆ in the solvent of choice, which should be purged with an inert gas. Rhein-¹³C₆ is soluble in DMSO.

Description

Rhein-¹³C₆ is intended for use as an internal standard for the quantification of rhein (Item No. 17345) by GC- or LC-MS. Rhein is an anthraquinone derivative that has been found in *R. rhubarbarum* and is an active metabolite of diacerein (Item No. 11710) that has diverse biological activities, including anticancer, antioxidant, and anti-inflammatory activities.¹⁻⁴ It induces cell cycle arrest at the S phase and inhibits the proliferation of HepG2 cells when used at concentrations of 40 and 100 μM, respectively.¹ Rhein (100 and 200 μM) also inhibits the proliferation of MCF-7 and MDA-MB-435s breast cancer cells under normoxic and hypoxic conditions.² It reduces controlled cortical impact-induced decreases in catalase and superoxide dismutase (SOD) activity and malondialdehyde (MDA), glutathione, and glutathione disulfide levels in the brain in a rat model of traumatic brain injury when administered at a dose of 12 mg/kg.³ Rhein (50 mg/kg per day) reduces cerulein-induced increases in serum TNF-α, IL-1β, and amylase levels, as well as reduces pancreatic glandular atrophy and fibrosis, in a mouse model of chronic pancreatitis.⁴

References

1. Liu, S., Wang, J., Shao, T., *et al.* The natural agent rhein induces β-catenin degradation and tumour growth arrest. *J. Cell. Mol. Med.* **22**(1), 589-599 (2018).
2. Fernand, V.E., Losso, J.N., Traux, R.E., *et al.* Rhein inhibits angiogenesis and the viability of hormone-dependent and -independent cancer cells under normoxic or hypoxic conditions *in vitro*. *Chem. Biol. Interact.* **192**(3), 220-232 (2011).
3. Xu, X., Lv, H., Xia, Z., *et al.* Rhein exhibits antioxidative effects similar to Rhubarb in a rat model of traumatic brain injury. *BMC Complement. Altern. Med.* **17**(1), 140 (2017).
4. Tsang, S.W., Zhang, H., Lin, C., *et al.* Rhein, a natural anthraquinone derivative, attenuates the activation of pancreatic stellate cells and ameliorates pancreatic fibrosis in mice with experimental chronic pancreatitis. *PLoS One* **8**(12), e82201 (2013).

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, 10/30/2024

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