

PRODUCT INFORMATION



Troxeutin

Item No. 27220

CAS Registry No.: 7085-55-4

Formal Name: 2-[3,4-bis(2-hydroxyethoxy)phenyl]-3-[[6-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranosyl]oxy]-5-hydroxy-7-(2-hydroxyethoxy)-4H-1-benzopyran-4-one

Synonyms:

MF: $C_{33}H_{42}O_{19}$

FW: 742.7

Purity: $\geq 90\%$

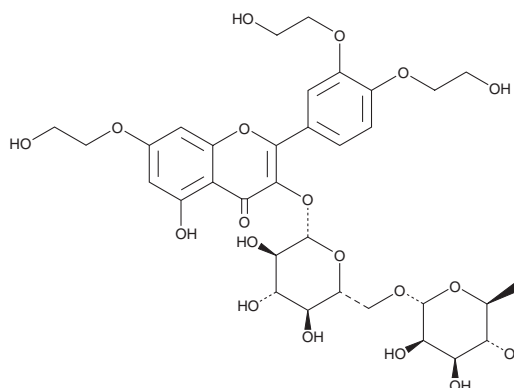
UV/Vis.: λ_{max} : 257, 352 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/Flos Sophorae Immaturus



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

Laboratory Procedures

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of troxeutin can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of troxeutin in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Troxeutin is a flavonoid that has been found in a variety of fruits and vegetables and has diverse biological activities.¹⁻³ It reduces D-galactose-induced increases in renal malondialdehyde (MDA) levels and decreases in renal Cu/Zn superoxide dismutase (SOD), catalase, and glutathione peroxidase (GPX) activities in mice when administered at a dose of 150 mg/kg per day, as well as attenuates D-galactose-induced increases in renal levels of NF- κ B p65, inducible nitric oxide synthase (iNOS), COX-2, and prostaglandin E₂ receptor 2 (EP₂).¹ Troxeutin (150 mg/kg per day) reduces blood glucose levels, plasma triglyceride levels, heart rate, and blood pressure in a rat model of type 2 diabetes induced by streptozotocin (STZ; Item No. 13104).² Troxeutin (150 mg/kg per day) prevents the loss of dopaminergic neurons in the substantia nigra pars compacta (SNc) and reduces apomorphine-induced contralateral rotations in a rat model of Parkinson's disease induced by 6-OHDA (Item No. 25330).³

References

1. Fan, S.-h., Zhang, Z.-f., Zheng, Y.-l., et al. *Int. Immunopharmacol.* **9**(1), 91-96 (2009).
2. Yu, Y. and Zheng, G. *Mol. Med. Rep.* **15**(6), 3473-3478 (2017).
3. Baluchnejadmojarad, T., Jamali-Raeufy, N., Zabihnejad, S., et al. *Eur. J. Pharmacol.* **801**, 72-78 (2017).

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, 11/11/2022

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