

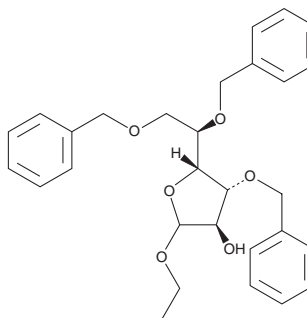
PRODUCT INFORMATION



Tribenoside

Item No. 33498

CAS Registry No.: 10310-32-4
Formal Name: ethyl 3,5,6-tris-O-(phenylmethyl)-D-glucofuranoside
MF: $C_{29}H_{34}O_6$
FW: 478.6
Purity: $\geq 95\%$ (mixture of anomers)
Supplied as: A neat oil
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Tribenoside is a vasoprotective agent with anti-allergic and anti-inflammatory activities.^{1,2} It inhibits agonist-induced contractions in various isolated porcine venous preparations when used at concentrations ranging from 10 to 30 $\mu\text{g/ml}$.¹ Tribenoside inhibits compound 48/80-induced histamine release in isolated rat peritoneal cells ($\text{EC}_{50} = 10 \mu\text{g/ml}$).³ It reduces decreases in proteoglycan levels and the severity of cartilage destruction in the inflamed joints in a rabbit model of surgically induced osteoarthritis when administered at doses of 100 and 200 mg/kg per day.⁴ Formulations containing tribenoside have been used in the treatment of hemorrhoids and varicose veins.

References

1. Jaques, R. The activity pattern, in isolated veins, of a benzyl glucofuranoside derivative possessing combined anti-inflammatory and venotropic effects. *Pharmacology* **4**(4), 193-202 (1970).
2. Jaques, R. The pharmacological activity of tribenoside. *Pharmacology* **15**(5), 445-460 (1977).
3. Rüegg, M. and Jaques, R. Tribenoside as an inhibitor of chemically induced histamine release. *Experientia* **30**(4), 399-401 (1974).
4. Colombo, C., Butler, M., Hickman, L., et al. A new model of osteoarthritis in rabbits. II. Evaluation of anti-osteoarthritic effects of selected antirheumatic drugs administered systemically. *Arthritis Rheum.* **26**(9), 1132-1139 (1983).

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, 12/07/2022

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