

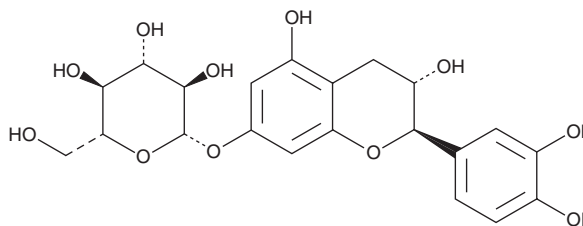
PRODUCT INFORMATION



Catechin 7-O-β-D-Glucopyranoside

Item No. 45760

CAS Registry No.: 65597-47-9
Formal Name: (2R,3S)-2-(3,4-dihydroxyphenyl)-3,4-dihydro-3,5-dihydroxy-2H-1-benzopyran-7-yl β-D-glucopyranoside
MF: C₂₁H₂₄O₁₁
FW: 452.4
Purity: ≥98%
Supplied as: A 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

Catechin 7-O-β-D-glucopyranoside is supplied as a solid. A stock solution may be made by dissolving the catechin 7-O-β-D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. Catechin 7-O-β-D-glucopyranoside is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Catechin 7-O-β-D-glucopyranoside is a flavonol glycoside that has been found in *P. longifolia* and has diverse biological activities.¹⁻³ It inhibits lipid peroxidation and scavenges hydroxyl radicals and superoxide anions in cell-free assays in a concentration-dependent manner and inhibits hydrogen peroxide-induced DNA damage and apoptosis in BJAB B cell lymphoma cells at 200 μM.¹ Catechin 7-O-β-D-glucopyranoside also inhibits lipid accumulation in 3T3-L1 cells.³ It decreases disease severity in a rat model of colitis at 2-10 mg/kg.³

References

1. Baek, J.A., Son, Y.-O., Fang, M., *et al.* Catechin-7-O-β-D-glucopyranoside scavenges free radicals and protects human B lymphoma BJAB cells on H₂O₂-mediated oxidative stress. *Food Sci. Biotechnol.* **20**(1), 151-158 (2011).
2. Wang, Z., Cui, Y., Xu, S., *et al.* Anti-lipid accumulation effects and mechanism of the extract and isolated compounds from *Potentilla longifolia*. *Fitoterapia* **168**, 105524 (2023).
3. Kook, S.-H., Choi, K.-C., Cho, S.-W., *et al.* Catechin-7-O-β-D-glucopyranoside isolated from the seed of *Phaseolus calcaratus* Roxburgh ameliorates experimental colitis in rat. *Int. Immunopharmacol.* **29**(2), 521-527 (2015).

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, 07/17/2026

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