

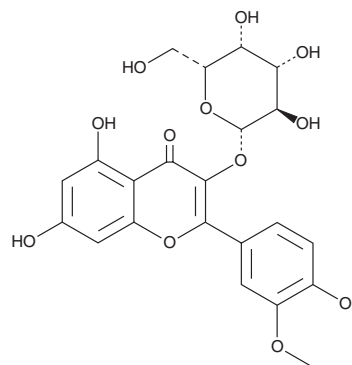
PRODUCT INFORMATION



Isorhamnetin-3-O-galactoside

Item No. 45821

CAS Registry No.: 6743-92-6
Formal Name: 3-(β-D-galactopyranosyloxy)-5,7-dihydroxy-2-(4-hydroxy-3-methoxyphenyl)-4H-1-benzopyran-4-one
Synonyms: Cacticin, I3G, IMG, Quercetin 3-methoxy-3-O-galactoside
MF: C₂₂H₂₂O₁₂
FW: 478.4
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

Isorhamnetin-3-O-galactoside is supplied as a solid. A stock solution may be made by dissolving the isorhamnetin-3-O-galactoside in the solvent of choice, which should be purged with an inert gas. Isorhamnetin-3-O-galactoside is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Isorhamnetin-3-O-galactoside is a flavonoid glycoside that has been found in *O. javanica* and has diverse biological activities.¹⁻³ It reduces thrombin and Factor Xa levels in human umbilical vein endothelial cells (HUVECs), increases activated partial thromboplastin time (aPTT) and prothrombin time in human plasma, and increases tail bleeding time in mice.² Isorhamnetin-3-O-galactoside (100 mg/kg) increases TNF-α serum levels and alanine transaminase (ALT) and aspartate aminotransferase (AST) serum activity and reduces malondialdehyde (MDA), inducible nitric oxide synthase (iNOS), and COX-2 serum levels in a mouse model of hepatic injury.³ It reduces peritoneal barrier permeability induced by high mobility group protein 1 (HMGB1) in mice and increases survival in a mouse model of sepsis.¹

References

1. Kim, T.H., Ku, S.-K., and Bae, J.-S. Anti-inflammatory activities of isorhamnetin-3-O-galactoside against HMGB1-induced inflammatory responses in both HUVECs and CLP-induced septic mice. *J. Cell. Biochem.* **114**(2), 336-345 (2013).
2. Ku, S.-K., Kim, T.H., Lee, S., et al. Antithrombotic and profibrinolytic activities of isorhamnetin-3-O-galactoside and hyperoside. *Food Chem. Toxicol.* **53**, 197-204 (2013).
3. Kim, D.-W., Cho, H.-I., Kim, K.-M., et al. Isorhamnetin-3-O-galactoside protects against CCl₄-induced hepatic injury in mice. *Biomol. Ther. (Seoul)* **20**(4), 406-412 (2012).

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, 08/07/2026

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