

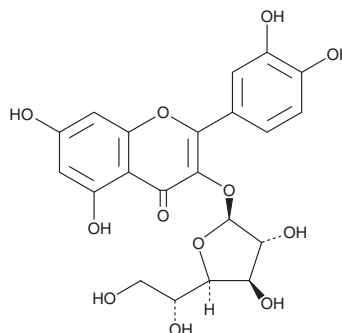
PRODUCT INFORMATION



Isoquercitroside

Item No. 20802

CAS Registry No.: 21637-25-2
Formal Name: 2-(3,4-dihydroxyphenyl)-3-(β-D-glucofuranosyloxy)-5,7-dihydroxy-4H-1-benzopyran-4-one
Synonyms: Quercetin 3-(β-D-glucofuranoside), Quercetin 3-O-β-D-glucofuranoside
MF: C₂₁H₂₀O₁₂
FW: 464.4
Purity: ≥98%
UV/Vis.: λ_{max}: 257, 361 nm
Supplied as: A crystalline 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

Isoquercitroside is supplied as a crystalline solid. A stock solution may be made by dissolving the isoquercitroside in the solvent of choice, which should be purged with an inert gas. Isoquercitroside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of isoquercitroside in these solvents is approximately 1, 30, and 15 mg/ml, respectively.

Isoquercitroside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, isoquercitroside should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Isoquercitroside has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Isoquercitroside is a 3-O-glucoside of quercetin (Item No. 10005169), that can be found in several plants. Isoquercitroside has higher bioavailability than quercetin and displays antioxidant and chemoprotective effects both *in vitro* and *in vivo*.¹

Reference

1. Valentová, K., Vrba, J., Baniřová, M., *et al.* Isoquercitrin: Pharmacology, toxicology, and metabolism. *Food Chem. Toxicol.* **68**, 267-282 (2014).

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, 12/01/2022

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