

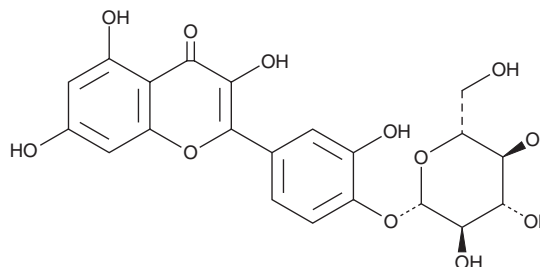
# PRODUCT INFORMATION



## Spiraeoside

Item No. 45667

**CAS Registry No.:** 20229-56-5  
**Formal Name:** 2-[4-(β-D-glucopyranosyloxy)-3-hydroxyphenyl]-3,5,7-trihydroxy-4H-1-benzopyran-4-one  
**Synonyms:** Quercetin-4-O-β-D-glycoside, Quercetin 4'-O-glucoside  
**MF:** C<sub>21</sub>H<sub>20</sub>O<sub>12</sub>  
**FW:** 464.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

Spiraeoside is supplied as a solid. A stock solution may be made by dissolving the spiraeoside in the solvent of choice, which should be purged with an inert gas. Spiraeoside is soluble (≥10 mg/ml) in ethanol and DMSO.

### Description

Spiraeoside is a flavonoid glycoside that has been found in *F. ulmaria* and has diverse biological activities.<sup>1,2</sup> It scavenges DPPH (Item No. 14805) and hydroxyl radicals (IC<sub>50</sub>s = 19.54 and 5.1 μg/ml, respectively) and inhibits β-glucuronidase and lipoxygenase (LO) in cell-free assays. Spiraeoside (15 μg/ml) induces cell cycle arrest and apoptosis in HeLa cervical cancer cells.<sup>2</sup> It decreases the severity of ethanol-induced gastric lesions in rats when administered at a dose of 50 mg/kg.<sup>1</sup>

### References

1. Samardžić, S., Arsenijević, J., Božić, D., *et al.* Antioxidant, anti-inflammatory and gastroprotective activity of *Filipendula ulmaria* (L.) Maxim. and *Filipendula vulgaris* Moench. *J. Ethnopharmacol.* **213**, 132-137 (2018).
2. Nile, A., Nile, S.H., Cespedes-Acuña, C.L., *et al.* Spiraeoside extracted from red onion skin ameliorates apoptosis and exerts potent antitumor, antioxidant and enzyme inhibitory effects. *Food Chem. Toxicol.* **154**, 112327 (2021).

#### 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, 05/22/2026

#### 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