

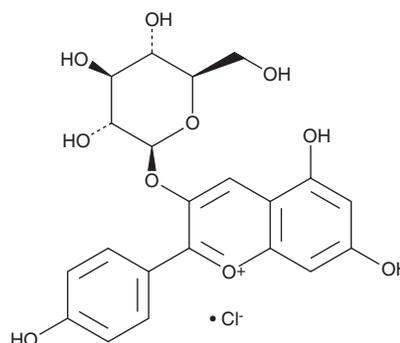
PRODUCT INFORMATION



Pelargonidin 3-O-glucoside (chloride)

Item No. 30224

CAS Registry No.: 18466-51-8
Formal Name: 3-(β-D-glucopyranosyloxy)-5,7-dihydroxy-2-(4-hydroxyphenyl)-1-benzopyrylium, monochloride
Synonym: Callistephin
MF: C₂₁H₂₁O₁₀ • Cl
FW: 468.8
Purity: ≥95%
UV/Vis.: λ_{max}: 272, 434, 517 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Vaccinium myrtillus*



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

Laboratory Procedures

Pelargonidin 3-O-glucoside (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the pelargonidin 3-O-glucoside (chloride) in the solvent of choice, which should be purged with an inert gas. Pelargonidin 3-O-glucoside (chloride) is soluble in chloroform, DMSO, ethyl acetate, and dichloromethane.

Description

Pelargonidin 3-O-glucoside is an anthocyanin and the 3-O-glucoside of pelargonidin (Item No. 19753) that has been found in *Fragaria* and has transporter inhibitory and antioxidant properties.^{1,2} It inhibits bilitranslocase transporter activity in isolated rat liver plasma membrane vesicles (K_i = 2.8 μM).¹ Pelargonidin 3-O-glucoside scavenges peroxy radicals in an oxygen radical absorbance capacity (ORAC) assay.²

References

1. Passamonti, S., Vrhovsek, U., and Mattivi, F. The interaction of anthocyanins with bilitranslocase. *Biochem. Biophys. Res. Commun.* **296**(3), 631-636 (2002).
2. Fukumoto, L.R. and Mazza, G. Assessing antioxidant and prooxidant activities of phenolic compounds. *J. Agric. Food. Chem.* **48**(8), 3597-3604 (2000).

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, 10/28/2022

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