

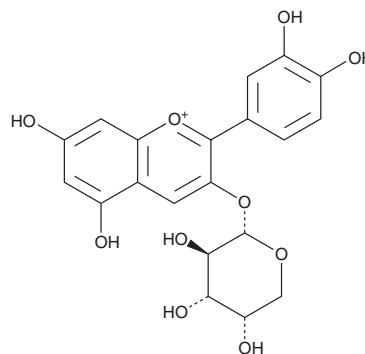
PRODUCT INFORMATION



Cyanidin 3-O-arabinoside

Item No. 26183

CAS Registry No.: 792868-19-0
Formal Name: 3-(α -L-arabinopyranosyloxy)-2-(3,4-dihydroxyphenyl)-5,7-dihydroxy-1-benzopyrylium
MF: C₂₀H₁₉O₁₀
FW: 419.4
Purity: \geq 95%
UV/Vis.: λ_{max} : 284, 539 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Cyanidin 3-O-arabinoside is supplied as a crystalline solid. A stock solution may be made by dissolving the cyanidin 3-O-arabinoside in the solvent of choice, which should be purged with an inert gas. Cyanidin 3-O-arabinoside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of cyanidin 3-O-arabinoside in these solvents is approximately 3.3, 10, and 17 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of cyanidin 3-O-arabinoside can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of cyanidin 3-O-arabinoside in PBS, pH 7.2, is approximately 0.14 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cyanidin 3-O-arabinoside is an anthocyanin antioxidant that has been found in eggplant.¹ It scavenges radicals in a Trolox equivalent antioxidant capacity (TEAC) assay.

Reference

1. Jing, P., Zhao, S., Ruan, S., *et al.* Quantitative studies on structure-ORAC relationships of anthocyanins from eggplant and radish using 3D-QSAR. *Food Chem.* **145**, 365-371 (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, 10/24/2022

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