

PRODUCT INFORMATION



Procyanidin

Item No. 29763

CAS Registry No.: 20347-71-1
Formal Name: 2-(3,4-dihydroxyphenyl)-2-[[2-(3,4-dihydroxyphenyl)-3,4-dihydro-5,7-dihydroxy-2H-1-benzopyran-3-yl]oxy]-3,4-dihydro-2H-1-benzopyran-3,4,5,7-tetrol

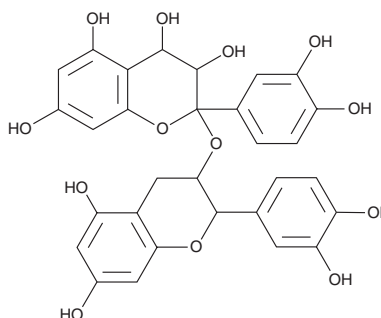
MF: C₃₀H₂₆O₁₃
FW: 594.5

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/Grape seeds



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

Laboratory Procedures

Procyanidin is supplied as a solid. A stock solution may be made by dissolving the procyanidin in the solvent of choice, which should be purged with an inert gas. Procyanidin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of procyanidin in these solvents is approximately 0.5, 10, and 3 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 procyanidin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of procyanidin in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Procyanidin is a polyphenol that exhibits antioxidant activity in a macrocyclic nickel complex-catalyzed Briggs-Rauscher oscillation system.¹

Reference

1. Li, M., Hu, G., and Chen, Y. Evaluation of the antioxidant capacity of natural polyphenolic compounds using a macrocyclic Ni-(II) complex-catalysed Briggs-Rauscher reaction. *Food Chem.* **197(Pt A)**, 987-991 (2015).

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, 10/06/2022

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