

PRODUCT INFORMATION

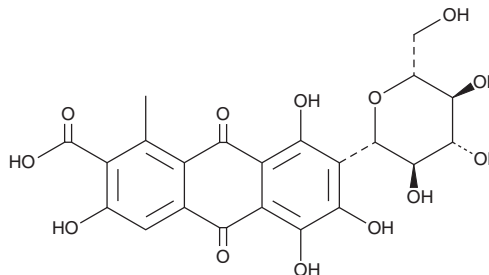


Cochineal

Item No. 35334

CAS Registry No.: 1260-17-9
Formal Name: 7-β-D-glucopyranosyl-9,10-dihydro-3,5,6,8-tetrahydroxy-1-methyl-9,10-dioxo-2-anthracenecarboxylic acid
Synonyms: Carminic Acid, Carmine 5297, E 120, NSC 326224

MF: C₂₂H₂₀O₁₃
FW: 492.4
Purity: ≥80%
UV/Vis.: λ_{max}: 223, 278 nm
Supplied as: A 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

Cochineal is supplied as a solid. A stock solution may be made by dissolving the cochineal in the solvent of choice, which should be purged with an inert gas. Cochineal is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

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 cochineal can be prepared by directly dissolving the solid in aqueous buffers. The solubility of cochineal in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cochineal is a hydroxyanthraquinone dye originally isolated from female cochineal insects (*D. coccus*).¹ Formulations containing cochineal have been used as color additives in food and pharmaceutical preparations.

Reference

1. Zuchelli, M., Villarruel, F.D., David-Gara, P., *et al.* Photophysics and photochemistry of carminic acid and related natural pigments. *Phys. Chem. Chem. Phys.* **22(17)**, 9534-9542 (2020).

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, 09/30/2022

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