

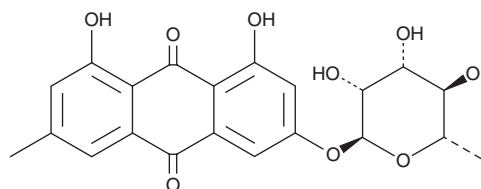
PRODUCT INFORMATION



Frangulin A

Item No. 36087

CAS Registry No.: 521-62-0
Formal Name: 3-[[6-deoxy- α -L-mannopyranosyl]oxy]-1,8-dihydroxy-6-methyl-9,10-anthracenedione
Synonym: Franguloside
MF: C₂₁H₂₀O₉
FW: 416.4
Purity: \geq 95%
UV/Vis.: λ_{max} : 225, 264, 280, 287, 423 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/Unknown Sp.



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

Laboratory Procedures

Frangulin A is supplied as a solid. A stock solution may be made by dissolving the frangulin A in the solvent of choice, which should be purged with an inert gas. Frangulin A is slightly soluble in ethanol and methanol.

Description

Frangulin A is an anthraquinone glycoside that has been found in *Rhamnus procumbens*.¹

Reference

1. Dwivedi, S.P.D., Pandey, V.B., Shah, A.H., *et al.* Chemical constituents of *Rhamnus procumbens* and pharmacological actions of emodin. *Phytother. Res.* **2(1)**, 51-53 (1988).

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, 12/19/2022

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