

PRODUCT INFORMATION



Sennidin A

Item No. 35237

CAS Registry No.: 641-12-3
Formal Name: (9R,9'R)-rel-(+)-9,9',10,10'-tetrahydro-4,4',5,5'-tetrahydroxy-10,10'-dioxo-[9,9'-bianthracene]-2,2'-dicarboxylic acid

MF: C₃₀H₁₈O₁₀

FW: 538.5

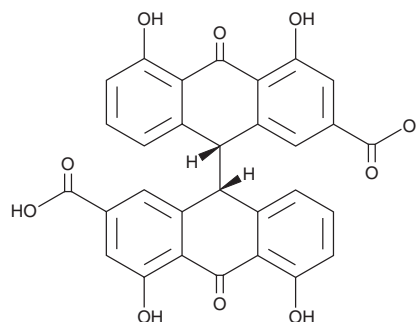
Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Plant/*Sennae folium*



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

Laboratory Procedures

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

Description

Sennidin A is an active metabolite of the dianthrone glycoside laxative sennoside A (Item No. 24969) that has also been found in *Senna*.^{1,2} It is formed from sennoside A by gut microbiota.¹ Sennidin A increases fecal excretion in mice in a dose-dependent manner.³

References

1. Hattori, M., Kim, G., Motoike, S., *et al.* Metabolism of sennosides by intestinal flora. *Chem. Pharm. Bull. (Tokyo)* **30(4)**, 1338-1346 (1982).
2. Metzger, W. and Reif, K. Determination of 1,8-dihydroxyanthranoids in senna. *J. Chromatogr. A* **740**, 133-138 (1996).
3. Fairbairn, J.W. and Moss, M.J. The relative purgative activities of 1,8-dihydroxyanthracene derivatives. *J. Pharm. Pharmacol.* **22(8)**, 584-593 (1970).

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, 10/05/2022

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