

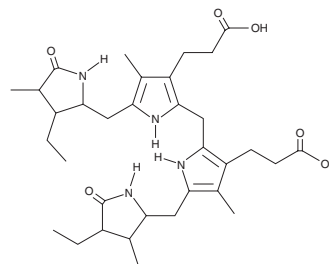
PRODUCT INFORMATION



Stercobilinogen (technical grade)

Item No. 36510

CAS Registry No.: 17095-63-5
Formal Name: (2R,3R,4S,16S,17R,18R)-2,17-diethyl-1,2,3,4,5,10,15,16,17,18,19,22,23,24-tetradecahydro-3,7,13,18-tetramethyl-1,19-dioxo-21H-bilene-8,12-dipropanoic acid
MF: C₃₃H₄₈N₄O₆
FW: 596.8
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

Stercobilinogen (technical grade) is supplied as a solid. A stock solution may be made by dissolving the stercobilinogen (technical grade) in the solvent of choice, which should be purged with an inert gas. Stercobilinogen (technical grade) is slightly soluble in methanol and DMSO.

Description

Stercobilinogen is a metabolite of bilirubin (Item No. 17161) and a precursor in the biosynthesis of the fecal pigment stercobilin (Item No. 36509).¹ Fecal levels of stercobilinogen are increased in patients with β -thalassemia.²

References

1. Vitek, L., Majer, F., Muchová, L., *et al.* Identification of bilirubin reduction products formed by *Clostridium fringens* isolated from human neonatal fecal flora. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **833(2)**, 149-157 (2006).
2. Gallo, E., Pich, P., Ricco, G., *et al.* The relationship between anemia, fecal stercobilinogen, erythrocyte survival, and globin synthesis in heterozygotes for β -thalassemia. *Blood* **46(5)**, 693-698 (1975).

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