

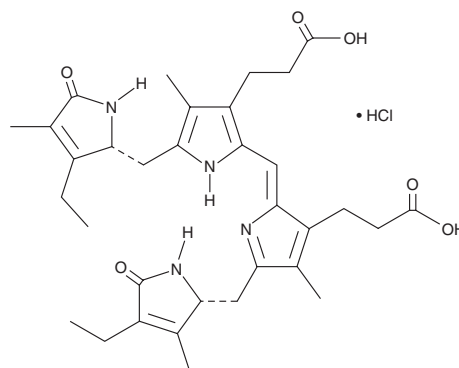
PRODUCT INFORMATION



Urobilin (hydrochloride)

Item No. 36508

CAS Registry No.: 28925-89-5
Formal Name: (4R,16R)-3,18-diethyl-1,4,5,15,16,19,22,24-octahydro-2,7,13,17-tetramethyl-1,19-dioxo-21H-bilane-8,12-dipropanoic acid, monohydrochloride
Synonym: Urochrome
MF: $C_{33}H_{42}N_4O_6 \cdot HCl$ (for d-Urobilin IX)
FW: 627.2
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

Urobilin (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the urobilin (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Urobilin (hydrochloride) is soluble in 1 M sodium hydroxide.

Description

Urobilin is a byproduct of hemoglobin degradation.¹ It is formed via oxidation of the bilirubin metabolite urobilinogen and is excreted in both urine and feces by mammals, including humans. Urobilin has been used as a marker of human waste contamination in recreational and source waters.

Reference

1. Jones-Lepp, T.L. Chemical markers of human waste contamination: Analysis of urobilin and pharmaceuticals in source waters. *J. Environ. Monit.* **8**(4), 472-478 (2006).

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, 04/03/2023

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