

PRODUCT INFORMATION



Biliverdin (hydrochloride)

Item No. 19257

CAS Registry No.: 856699-18-8
Formal Name: 3,18-diethenyl-1,19,22,24-tetrahydro-2,7,13,17-tetramethyl-1,19-dioxo-21H-biliverdin-8,12-dipropanoic acid, monohydrochloride

MF: C₃₃H₃₄N₄O₆ • HCl
FW: 619.1

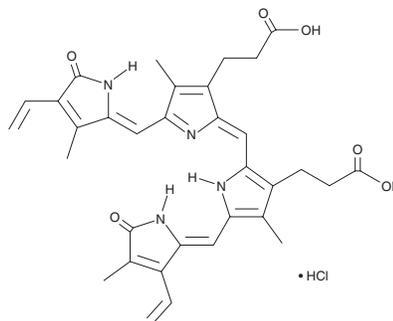
Purity: ≥95%

UV/Vis.: λ_{max}: 310, 376, 689 nm

Supplied as: A crystalline 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

Biliverdin (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the biliverdin (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Biliverdin (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of biliverdin (hydrochloride) in these solvents is approximately 20 mg/ml.

Biliverdin (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, biliverdin (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Biliverdin (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Biliverdin is a green bile pigment produced from the oxidation of heme in a reaction catalyzed by heme oxygenase and is further reduced to bilirubin (Item No. 17161) by biliverdin reductase.^{1,2} Biliverdin regulates the cellular heme degradation process by inhibiting substrates from binding to the catalytic site of heme oxygenase. Bile pigments such as biliverdin are known to possess anti-mutagenic and antioxidant properties^{3,4}

References

1. Liu, Y., Moënné-Loccoz, P., Loehr, T.M., *et al.* Heme oxygenase-1, intermediates in verdoheme formation and the requirement for reduction equivalents. *J. Biol. Chem.* **272**, 6909-6917 (1997).
2. Maines, M.D. The heme oxygenase system: A regulator of second messenger gases. *Annu. Rev. Pharmacol. Toxicol.* **37**, 517-554 (1997).
3. Bulmer, A.C., Ried, K., Blanchfield, J.T., *et al.* The anti-mutagenic properties of bile pigments. *Mutat. Res.* **658(1-2)**, 28-41 (2008).
4. Mölzer, C., Huber, H., Steyrer, A., *et al.* *In vitro* antioxidant capacity and antigenotoxic properties of protoporphyrin and structurally related tetrapyrroles. *Free Rad. Res.* **46(11)**, 1369-1377 (2012).

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, 03/09/2023

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