

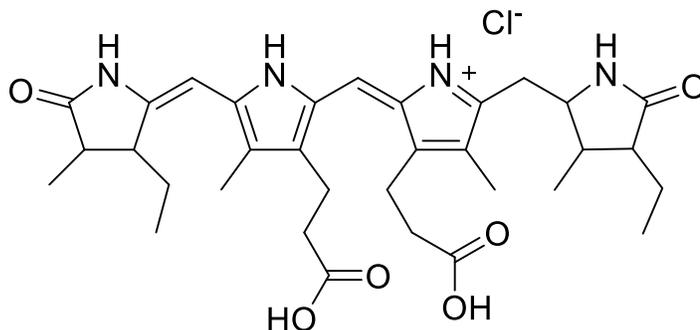
Stercobilin Hydrochloride

Catalog number: F-H150

Molecular Formula: $C_{33}H_{47}N_4O_6Cl$

MW: 631.22

CAS: 34217-90-8



Solubility: DMF, pyridine and basic solutions > pH 9.0

Storage and Handling: Stercobilin Hydrochloride is stable for at least one year when stored as a solid, protected from moisture, at $-20^{\circ}C$. Protect from light.

Background: Stercobilin Hydrochloride is produced via the oxidation of stercobilinogen, a degradation product of bilirubin, and excreted in the feces. Bilirubin is a water insoluble tetrapyrrole produced from the reduction of biliverdin in a reaction catalyzed by the enzyme bilirubin reductase. Water insoluble bilirubin (also called indirect bilirubin) in vivo undergoes glucuronidation in the liver (addition of one or two glucuronic acids through a glycosidic bond) to form the water soluble bilirubin mono or diglucuronide (also called bilirubin conjugate or direct bilirubin). Bilirubin conjugate is excreted from the liver in bile or is converted to mesobilinogen via gut bacteria and then to urobilinogen and excreted in the urine as urobilin or stercobilinogen and excreted in the feces as stercobilin. Stercobilin Hydrochloride is soluble in basic aqueous solutions (pH > 9 for initial dissolution) and soluble down to pH 7 once in solution as well as methanol and ethanol if made slightly basic.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

