

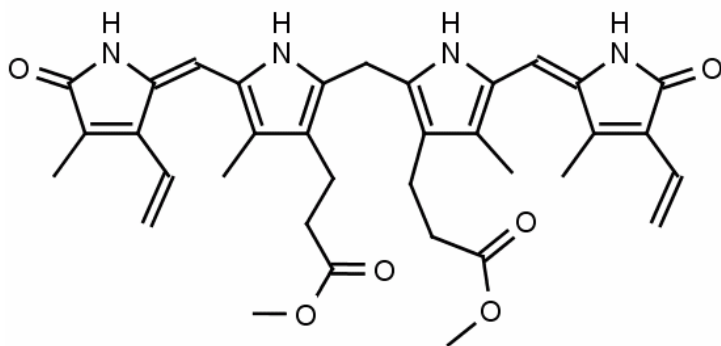


Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Product Name: Bilirubin dimethyl ester



Catalog Number	CAS Registry No.	MW (g/mol)	Chemical Formula
F-H110	19792-68-8	612.73	C ₃₅ H ₄₀ N ₄ O ₆

Storage: Bilirubin dimethyl ester is stable for at least one year when stored as a solid, protected from moisture, at -20 °C. Protect from light.

Field of Use: Bilirubin dimethyl ester is a natural derivative of bilirubin and is found in normal sera representing an average of 1.75% total sera bilirubin¹. 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 urobilinogen and excreted in the urine as urobilin or in the feces as stercobilin. Bilirubin dimethyl ester has been found to be converted to bilirubin conjugate via esterase and glucuronidase activity *in vivo*².

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.

References:

1. Muraca, M.; Blanckaert, N. Liquid-chromatographic assay and identification of mono- and diester conjugates of bilirubin in normal serum. *Clin Chem* **1983**, 29, 1767-71.
2. Burchell, B.; Odell, G. B. A rat liver microsomal carboxyesterase and a bilirubin UDP-glucuronyl transferase are responsible for the formation of bilirubin glucuronides from bilirubin dimethyl ester. *FEBS Lett* **1981**, 135, 304-8.

Echelon Biosciences products are sold for research and development purposes only and are not for diagnostic use or to be incorporated into products for resale without written permission from Echelon Biosciences. Materials in this publication, as well as applications and methods and use, may be covered by one or more U.S. or foreign patents or patents pending. We welcome inquiries about licensing the use of our trademarks and technologies at busdev@echelon-inc.com.