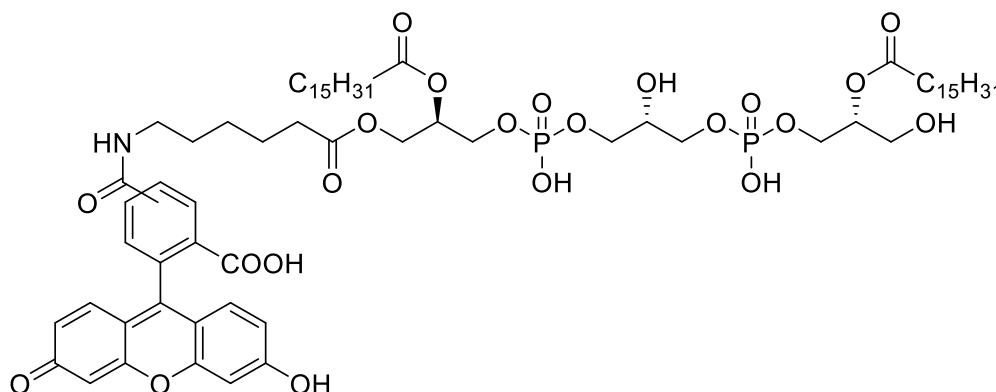


5(6)-Carboxyfluorescein-Labeled Monolysocardiolipin (Fluorescein-MCL)

Catalog number: L-M16F



Molecular Formula: C₆₈H₁₀₃NO₂₂P₂

MW: 1348.49

CAS: n/a

Solubility: Chloroform, methanol, ethanol >1mg/mL

Storage and Handling: Fluorescein-MCL is stable for at least one year when stored as a solid, protected from moisture, and light at -20 °C. Reconstitute with methanol and dilute with water or neutral pH, buffered salt solutions, i.e. PBS, TBS, etc. Should be stored in glass containers to prevent material loss due to adsorption to the vessel surface. Storage in basic (pH > 9) or acidic (pH < 4) buffers will result in slow decomposition of the product. After reconstitution, solutions of biotin-MCL should be stored at -20 °C between uses.

Background: Cardiolipin (CL) is an important component of the prokaryotic and eukaryotic inner mitochondrial membrane. In humans, CL is made by cardiolipin synthase from phosphatidylglycerol and cytidinediphosphate-diacylglycerol. Monolysocardiolipin is an intermediate in CL remodeling. Acetyl transferase tafazzin mutations impairing linoleic acid transfer to MCL are considered the main reason of Barth syndrome; and autoantibodies to MCL have been found in patients with anti-phospholipid syndrome and lupus. Biotinylated lipids have been used in microplate and bead-based assays and affinity experiments.

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.

Technical Data Sheet, Rev 2, 2-12-19 – For research use only. Not intended for diagnostic or therapeutic use.

