Leukotriene A4 methyl ester

Substrate for LTA_4 hydrolase and LTC_4 synthase

Unstable intermediate in the biosynthesis of LTB_4 and LTC_4 . The naturally occuring free acid is a substrate for LTA_4 hydrolase and LTC_4 synthase and plays a central role in transcellular leukotrieneand lipoxin biosynthesis. Mobilizes Ca^{2+} in human neutrophils. Supplied as the methyl ester for greater stability.

Citations: 4

View Online »

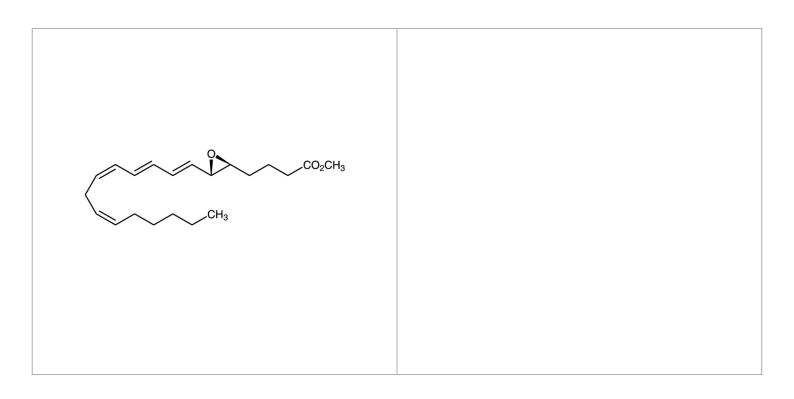
Ordering Information

Order Online »

BML-LA004-1000	1mg
BML-LA004-0050	50µg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored, as supplied, at -80°C.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name LTA4-ME

Appearance Liquid.

CAS 73466-12-3

Couple Target Leukotriene hydrolase, Leukotriene synthase, Prostanoid

receptor

Couple Type Ligand, Substrate

Formula $C_{21}H_{32}O_3$

Formulation Oil dissolved in 2% triethylamine/hexane.

MW 332.5

Purity ≥95% (HPLC) (remainder 11-trans isomer)

Source Synthetic.

Last modified: May 29, 2024

