8(S)-Hydroxyeicosa-5Z,9E,11Z,14Ztetraenoic acid (HETE)

Activator of PKC and PPARa

Biosynthesis in human neutrophils, phorbol ester-treated mouse epidermis and by hepatic microsomal cytochrome P-450. Chemotactic for neutrophils. Activator of PPARα. Activates mouse skin protein kinase C (PKC).

Citations: 4

View Online »

Ordering Information

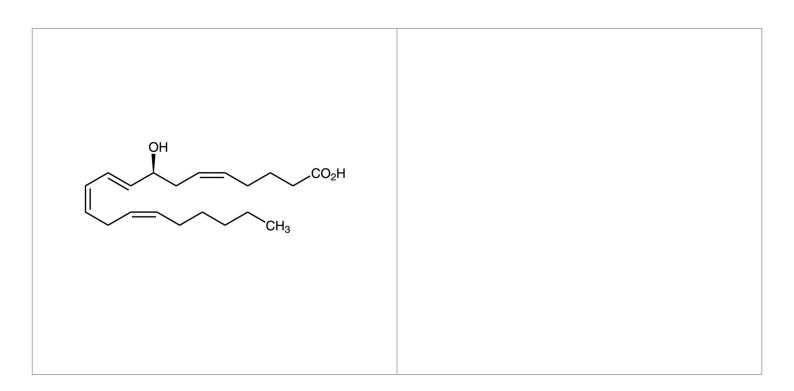
Order Online »

BML-H008-0050

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 8(S)-HETE

CAS 98462-03-4

Couple Target PKC, PPAR

Couple Type Activator

Formula $C_{20}H_{32}O_3$

Formulation Liquid. In ethanol.

MW 320.5

Purity ≥99% (HPLC)

Technical Info / Product NotesOpen ampule and aliquot into suitable tubes or vials. Store

at -80°C or colder. For experiments with cells or tissues, aliquots of ethanol solutions may be added directly to aqueous media. For HPLC, aliquots of ethanol solutions should be evaporated to dryness under a stream of nitrogen and HPLC mobile phase added to the desired

concentration.

Last modified: May 29, 2024

