

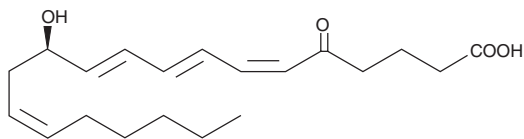
PRODUCT INFORMATION



5-oxo Leukotriene B₄

Item No. 35416

CAS Registry No.: 195061-92-8
Formal Name: 12R-hydroxy-5-oxo-6Z,8E,10E,14Z-eicosatetraenoic acid
Synonyms: 5-keto LTB₄, 5-oxo LTB₄
MF: C₂₀H₃₀O₄
FW: 334.5
Purity: ≥90%
UV/Vis.: λ_{max}: 313 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

5-oxo Leukotriene B₄ (5-oxo LTB₄) is a byproduct in the synthesis of 12-oxo LTB₄ (Item No. 20140).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/14/2021

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM