

PRODUCT INFORMATION



Cysteinyl Leukotriene HPLC Mixture II

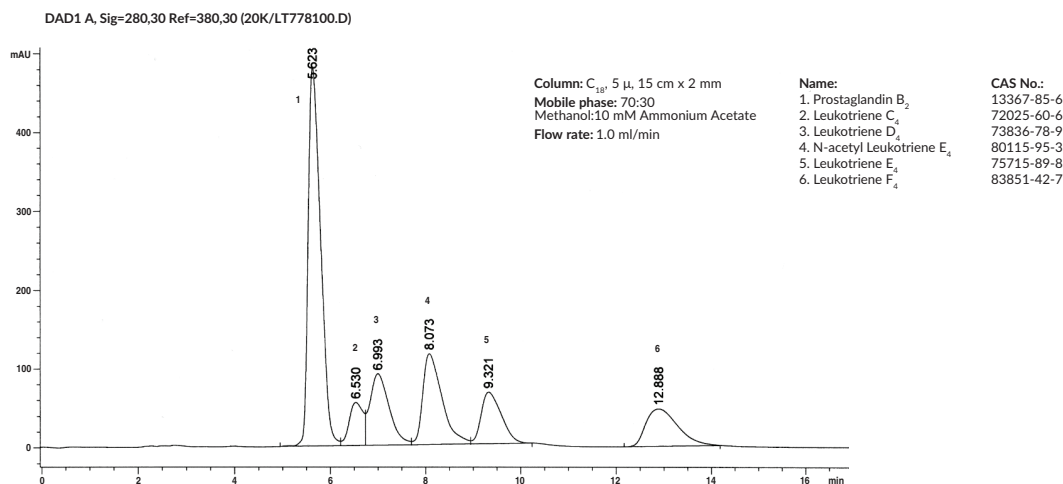
Item No. 20002

Purity: ≥98% for each compound
Supplied as: A solution in ethanol (Please see below for specific amounts of each compound)
Storage: -80°C
Stability: ≥6 months
Special Conditions: Light sensitive

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

Description and Data

This mixture contains all primary cysteinyl-leukotrienes derived from arachidonic acid and also includes Prostaglandin B₂. Contents: Leukotriene C₄, Leukotriene D₄, Leukotriene E₄, Leukotriene F₄, N-acetyl Leukotriene E₄ (5 µg each) and Prostaglandin B₂ (10 µg).



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, 08/04/2022

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