

# PRODUCT INFORMATION



## Vasoactive Eicosanoid HPLC Mixture

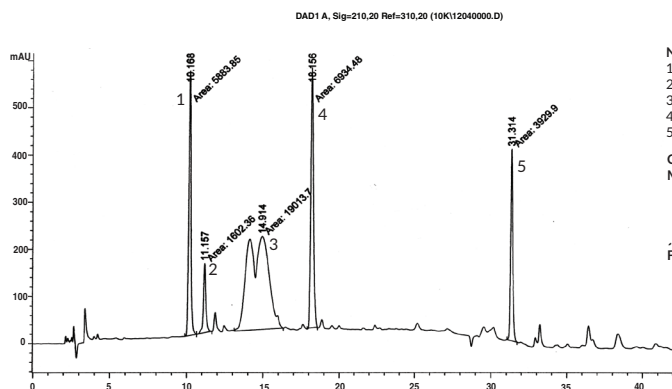
Item No. 10003

**Purity:** ≥98% for each compound  
**Supplied as:** A solution in methyl acetate (Please see below for specific amounts of each compound.)  
**Storage:** -20°C  
**Stability:** ≥6 months

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

### Laboratory Procedures

The vasoactive eicosanoid HPLC mixture contains a combination of Prostaglandin I<sub>2</sub> (PGI<sub>2</sub>) and Thromboxane A<sub>2</sub> (TXA<sub>2</sub>) supplied in methyl acetate. This HPLC mixture contains 5 µg of 12(S)-HHTrE and 100 µg of TXB<sub>2</sub>, 11-dehydro TXB<sub>2</sub>, 6-keto PGF<sub>1α</sub>, and 2,3-dinor-6-keto PGF<sub>1α</sub>.



**Name:**  
1. 6-keto Prostaglandin F<sub>1α</sub>  
2. 2,3-dinor-6-keto Prostaglandin F<sub>1α</sub>  
3. Thromboxane B<sub>2</sub>  
4. 11-dehydro Thromboxane B<sub>2</sub>  
5. 12(S)-HHTrE

**CAS No.:**  
58962-34-8  
54397-85-2  
67910-12-7  
54397-84-1

**Column:** C18, 5µm, 4.6 x 250 mm  
**Mobile phase:** Acetonitrile:Water:Trifluoroacetic Acid  
0 min. 25:75:0.1  
20 min 45:55:0.1  
30 min 65:35:0.1  
40 min 65:35:0.1

**Flow rate:** 1.0 ml/min

#### 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, 08/03/2022

#### 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