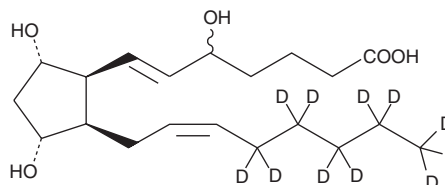


# PRODUCT INFORMATION



(±)5-iPF<sub>2α</sub>-VI-d<sub>11</sub>  
Item No. 10006654

**CAS Registry No.:** 936565-17-2  
**Formal Name:** (±)5,9α,11α-trihydroxy-(8β)-prosta-6E,14Z-dien-1-oic-16,16,17,17,18,18,19,19,20,20,20-d<sub>11</sub> acid  
**MF:** C<sub>20</sub>H<sub>23</sub>D<sub>11</sub>O<sub>5</sub>  
**FW:** 365.6  
**Chemical Purity:** ≥90% ((±)5-iPF<sub>2α</sub>-VI)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>11</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year



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

## Laboratory Procedures

(±)5-iPF<sub>2α</sub>-VI-d<sub>11</sub> contains 11 deuterium atoms at the 16, 16', 17, 17', 18, 18', 19, 19', 20, 20, and 20 positions. (±)5-iPF<sub>2α</sub>-VI-d<sub>11</sub> is intended for use as an internal standard for quantification of (±)5-iPF<sub>2α</sub>-VI (Item No. 16300) by GC- or LC-mass spectrometry (MS).

(±)5-iPF<sub>2α</sub>-VI-d<sub>11</sub> is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of (±)5-iPF<sub>2α</sub>-VI-d<sub>11</sub> in these solvents is approximately 50 mg/ml.

(±)5-iPF<sub>2α</sub>-VI-d<sub>11</sub> is used as an internal standard for the quantification of (±)5-iPF<sub>2α</sub>-VI by stable isotope dilution MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the weight indicated on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard (±)5-iPF<sub>2α</sub>-VI by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

## Description

Isoprostanes are prostaglandin (PG)-like products of free-radical induced lipid peroxidation.<sup>1</sup> Although the isoprostanes derived from arachidonic acid are the best characterized, many other polyunsaturated fatty acids can form isoprostanes.<sup>2</sup> iPF<sub>2α</sub>-VI is one of dozens of possible stereo- and regioisomeric isoprostanes which can be formed from arachidonic acid. To date, the most extensively studied of these is 8-isoprostane (8-*epi*-PGF<sub>2α</sub>; iPF<sub>2α</sub>-III).<sup>3,4</sup> However, 8-isoprostane is a minor isoprostane constituent when compared to some of the other isomers which form in natural conditions of oxidative stress,<sup>5</sup> including iPF<sub>2α</sub>-VI of the type-VI isoprostanes. This class has been shown to be one of the major isoprostane products, in contrast to 8-isoprostane. In addition to being produced in greater abundance than 8-isoprostane, type VI isoprostanes form internal lactones which facilitate their extraction and purification from biological samples.<sup>5-8</sup>

## References

1. Morrow, J.D., Hill, K.E., Burk, R.F., *et al. Proc. Natl. Acad. Sci. USA* **87**, 9383-9387 (1990).
2. Parchmann, S. and Mueller, M.J. *J. Biol. Chem.* **273**, 32650-32655 (1998).
3. Delanty, N., Reilly, M., Pratico, D., *et al. Br. J. Clin. Pharmacol.* **42**, 15-19 (1996).
4. Reilly, M.P., Barry, P., Lawson, J.A., *et al. Fibrinolysis & Proteolysis* **11**, 81-84 (1997).
5. Reilly, M.P., Pratico, D., Delanty, N., *et al. Circulation* **98**, 2822-2828 (1998).
6. Li, H., Lawson, J.A., Reilly, M., *et al. Proc. Natl. Acad. Sci. USA* **96**(23), 13381-13386 (1999).
7. Lawson, J.A., Li, H., Rokach, J., *et al. J. Biol. Chem.* **273**, 29295-29301 (1998).
8. Praticò, D., Barry, O.P., Lawson, J.A., *et al. Proc. Natl. Acad. Sci. USA* **95**, 3449-3454 (1998).

### 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, 07/13/2018

## 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