

# PRODUCT INFORMATION

## Butaprost (free acid)

Item No. 13741

**CAS Registry No.:** 433219-55-7

**Formal Name:** 9-oxo-11 $\alpha$ ,16S-dihydroxy-17-cyclobutyl-prost-13E-en-1-oic acid

**Synonyms:** ( $\pm$ )-15-deoxy-16S-hydroxy-17-cyclobutyl PGE<sub>1</sub>, 15-deoxy-16S-hydroxy-17-cyclobutyl PGE<sub>1</sub>

**MF:** C<sub>23</sub>H<sub>38</sub>O<sub>5</sub>

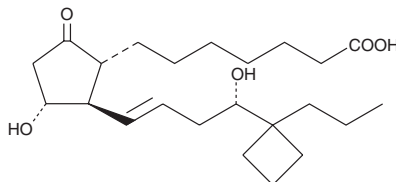
**FW:** 394.6

**Purity:**  $\geq 95\%$

**Supplied as:** A solution in methyl acetate

**Storage:** -20°C

**Stability:**  $\geq 2$  years



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

### Laboratory Procedures

Butaprost (free acid) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of butaprost (free acid) in ethanol is approximately 50 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Butaprost (free acid) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of butaprost (free acid) should be diluted with the aqueous buffer of choice. The solubility of butaprost (free acid) in PBS (pH 7.2) is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Butaprost is an EP<sub>2</sub> selective agonist which has frequently been used to pharmacologically define the EP receptor expression profile of various human and animal tissues and cells.<sup>1,2</sup> Prostaglandin free acids generally bind to their cognate receptors with 10 to 100 times the affinity of the corresponding ester derivative. Consistent with this trend, butaprost binds to membranes from EP<sub>2</sub> receptor-transfected CHO cells with a K<sub>i</sub> value of 2,400 nM, whereas butaprost (free acid) and CAY10399 (the 2-series congener of butaprost free acid) exhibit significantly lower K<sub>i</sub> values of 73 and 92 nM, respectively.<sup>3</sup> Butaprost (free acid) is therefore another useful tool for characterizing EP receptor-mediated signaling events.

### References

1. Kiriya, M., Ushikubi, F., Kobayashi, T., *et al.* Ligand binding specificities of the eight types and subtypes of the mouse prostanoid receptors expressed in Chinese hamster ovary cells. *Br. J. Pharmacol.* **122**(2), 217-224 (1997).
2. Lawrence, R.A. and Jones, R.L. Investigation of the prostaglandin E (EP-) receptor subtype mediating relaxation of the rabbit jugular vein. *Br. J. Pharmacol.* **105**(4), 817-824 (1992).
3. Tani, K., Naganawa, A., Ishida, A., *et al.* Design and synthesis of a highly selective EP<sub>2</sub>-receptor agonist. *Bioorg. Med. Chem. Lett.* **11**(15), 2025-2028 (2001).

#### 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, 06/29/2023

#### 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