

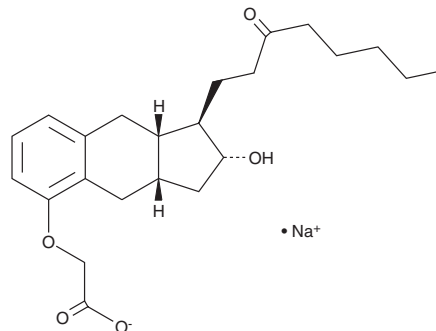
# PRODUCT INFORMATION



## 15-keto Treprostinil (sodium salt)

Item No. 22247

**Formal Name:** 2-(((1R,2R,3aS,9aS)-2-hydroxy-1-(3-oxooctyl)-2,3,3a,4,9,9a-hexahydro-1H-cyclopenta[b]naphthalen-5-yl)oxy)acetate, monosodium salt  
**MF:** C<sub>23</sub>H<sub>31</sub>O<sub>5</sub> • Na  
**FW:** 410.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 279 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

15-keto Treprostinil (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 15-keto treprostinil (sodium salt) in the solvent of choice. 15-keto Treprostinil (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 15-keto treprostinil (sodium salt) in ethanol and DMSO is approximately 10 mg/ml and approximately 20 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 15-keto treprostinil (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 15-keto treprostinil (sodium salt) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

15-keto Treprostinil is an impurity found in treprostinil (Item No. 10162), which is a stable analog of prostaglandin I<sub>2</sub> (prostacyclin; Item No. 18220) with a longer plasma half-life.<sup>1</sup> Formulations containing treprostinil are used for the treatment of primary pulmonary hypertension.

### Reference

1. Olschewski, H., Rose, F., Schermuly, R., *et al.* Prostacyclin and its analogues in the treatment of pulmonary hypertension. *Pharmacol. Ther.* **102**, 139-153 (2004).

#### 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, 06/09/2017

#### 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