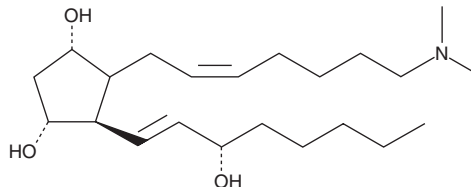


# PRODUCT INFORMATION



## Prostaglandin F<sub>2α</sub> dimethyl amine Item No. 16033

**CAS Registry No.:** 67508-09-2  
**Formal Name:** 1-(dimethylamino)-prosta-5Z,13E-diene-9α,11α,15S-triol  
**Synonyms:** Dinoprost dimethyl amine, PGF<sub>2α</sub> dimethyl amine  
**MF:** C<sub>22</sub>H<sub>41</sub>NO<sub>3</sub>  
**FW:** 367.6  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Prostaglandin F<sub>2α</sub> dimethyl amine (PGF<sub>2α</sub> dimethyl amine) 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 methyl acetate, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of PGF<sub>2α</sub> dimethyl amine in methyl acetate is approximately 10 mg/ml and approximately 50 mg/ml in DMSO and DMF.

PGF<sub>2α</sub> dimethyl amine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of PGF<sub>2α</sub> dimethyl amine should be diluted with the aqueous buffer of choice. The solubility of PGF<sub>2α</sub> dimethyl amine in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PGF<sub>2α</sub> dimethyl amine is a derivative of PGF<sub>2α</sub> which was designed as a PG antagonist for *in vitro* and *in vivo* studies.<sup>1</sup> PGF<sub>2α</sub> dimethyl amine is a weak FP receptor antagonist.<sup>1</sup> In gerbil colon, PGF<sub>2α</sub> dimethyl amine at a dose of 3.2 μg/ml inhibits the contractile effects of PGF<sub>2α</sub> (at 6 ng/ml) by 60%.<sup>1</sup>

### References

1. Maddox, Y.T., Ramwell, P.T., Shiner, C.S., *et al.* Amide and l-amino derivatives of F prostaglandins as prostaglandin antagonists. *Nature* **273**(5663), 549-552 (1978).

#### 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, 02/05/2024

#### 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