

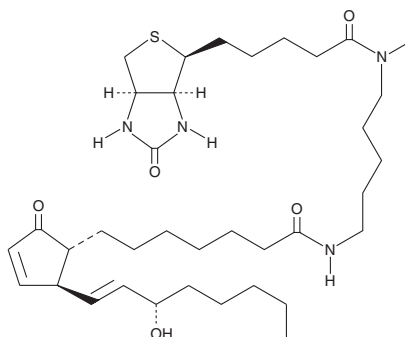
# PRODUCT INFORMATION



## Prostaglandin A<sub>1</sub>-biotin

Item No. 10013

**Formal Name:** N-9-oxo-15S-hydroxy-prosta-10,13E-dien-1-oyl-N'-biotinoyl-1,6-diaminopentane  
**Synonym:** PGA<sub>1</sub>-biotin  
**MF:** C<sub>35</sub>H<sub>58</sub>N<sub>4</sub>O<sub>5</sub>S  
**FW:** 646.9  
**Purity:** ≥95%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in ethanol



### Laboratory Procedures

For long term storage, we suggest that prostaglandin A<sub>1</sub>-biotin (PGA<sub>1</sub>-biotin) be stored as supplied at -20°C. It should be stable for at least one year.

PGA<sub>1</sub>-biotin 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of PGA<sub>1</sub>-biotin in these solvents is approximately 50, 25, and 50 mg/ml, respectively.

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. If an organic solvent-free solution of PGA<sub>1</sub>-biotin is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of PGA<sub>1</sub>-biotin in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

PGA<sub>1</sub> is one of the cyclopentenone prostaglandins, which have well documented antimitotic and antiproliferative effects.<sup>1,2</sup> The activity of the compounds in this class, which includes prostaglandins in both the A- and J-series, may result from changes in gene expression and the interaction with non-classical (*i.e.*, non-G protein-coupled receptor) pathways. PGA<sub>1</sub>-biotin is an affinity probe which allows PGA<sub>1</sub> to be detected through an interaction with the biotin ligand. PGA<sub>1</sub>-biotin was designed to allow PGA<sub>1</sub> to be detected in complexes with nucleic acid or protein binding partners. It is thus a tool to be used in the general elucidation of the mechanism of action of the cyclopentenone prostaglandins.

### References

1. Bregman, M.D. and Meyskens, F.L., Jr. Inhibition of human malignant melanoma colony-forming cells *in vitro* by prostaglandin A<sub>1</sub>. *Cancer Res.* **43**, 1642-1645 (1983).
2. Lical, P.M., Amici, C., Bonmassar, E., *et al.* Effects of cyclopentenone prostaglandins on myeloid cells during early infection with HTLV-I. II. Regulation of synthesis of inducible p72 heat shock protein. *J. Pharmacol. Exp. Ther.* **271**, 1096-1102 (1994).

#### 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, 07/29/2015

#### 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