

# PRODUCT INFORMATION



## Prostaglandin E<sub>2</sub> serinol amide

Item No. 10193

CAS Registry No.: 951209-67-9

Formal Name: N-[(2-hydroxy-1-hydroxymethyl)-ethyl]-9-oxo-11 $\alpha$ ,15S-dihydroxy-prosta-5Z,13E-dien-1-amide

Synonym: PGE<sub>2</sub>-SA

MF: C<sub>23</sub>H<sub>39</sub>NO<sub>6</sub>

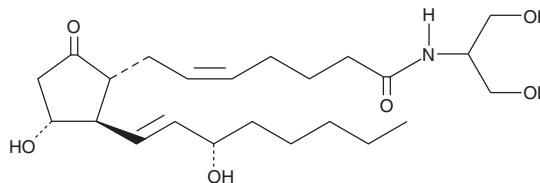
FW: 425.6

Purity:  $\geq$ 95%

Supplied as: A solution in ethanol

Storage: -80°C

Stability:  $\geq$ 2 years



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

### Laboratory Procedures

PGE<sub>2</sub>-SA 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 PGE<sub>2</sub>-SA in these solvents is approximately 30 mg/ml

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 ethanol is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of PGE<sub>2</sub>-SA in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

2-Arachidonyl glycerol (2-AG) exhibits cannabinoid agonist activity at the CB<sub>1</sub> receptor, is an important endogenous monoglyceride species, and is thus considered to be the natural ligand for the CB<sub>1</sub> receptor.<sup>1,2</sup> 2-AG can also be metabolized sequentially by COX-2 and specific PG synthases to form PG 2-glycerol esters.<sup>3</sup> PGE<sub>2</sub>-SA is a stable analog of PGE<sub>2</sub> 2-glycerol ester. The biological activity of PGE<sub>2</sub>-SA has not yet been determined.

### References

1. Sugiura, T., Kodaka, T., Kondo, S., *et al.* Is the cannabinoid CB<sub>1</sub> receptor a 2-arachidonoylglycerol receptor? Structural requirements for triggering a Ca<sup>2+</sup> transient in NG108-15 cells. *J. Biochem.* **122**, 890-895 (1997).
2. Kondo, S., Kondo, H., Nakane, S., *et al.* 2-Arachidonoylglycerol, and endogenous cannabinoid receptor agonist: Identification as one of the major species of monoacylglycerols in various rat tissues, and evidence for its generation through Ca<sup>2+</sup> -dependent and -independent mechanisms. *FEBS Lett.* **429**, 152-156 (1998).
3. Kozak, K.R., Crews, B.C., Morrow, J.D., *et al.* Metabolism of the endocannabinoids, 2-arachidonoylglycerol and anandamide, into prostaglandin, thromboxane, and prostacyclin glycerol esters and ethanolamides. *J. Biol. Chem.* **277**(47), 44877-44885 (2002).

#### 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, 03/07/2024

#### 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