

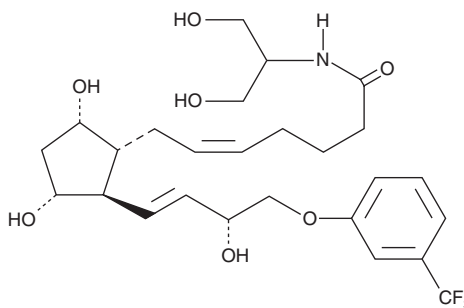
PRODUCT INFORMATION



Fluprostenol serinol amide

Item No. 10004236

CAS Registry No.: 1176658-85-7
Formal Name: N-[(2-hydroxy-1-hydroxymethyl)ethyl]-9 α ,11 α ,15R-trihydroxy-16-(3-trifluoromethyl)phenoxy)-17,18,19,20-tetranor-prosta-5Z,13E-dien-1-amide
Synonym: Flu-SA
MF: C₂₆H₃₆F₃NO₇
FW: 531.6
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 222, 277 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥ 1 year



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

Laboratory Procedures

Fluprostenol serinol amide 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 fluprostenol serinol amide in these solvents is approximately 20 and 30 mg/ml, respectively.

Fluprostenol serinol amide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of fluprostenol serinol amide should be diluted with the aqueous buffer of choice. The solubility of fluprostenol serinol amide 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 (CB) agonist activity at the CB₁ receptor,¹ is an important endogenous monoglyceride species,² and is thus considered to be the natural ligand for the CB₁ receptor. 2-AG can also be metabolized by cyclooxygenase-2 and specific prostaglandin H₂ (PGH₂) isomerases to form PG 2-glyceryl esters.³ Fluprostenol serinol amide (Flu-SA) is a stable analog of PGF_{2 α} 2-glyceryl ester that has much greater stability. The biological activity of Flu-SA has not yet been determined.

References

1. Sugiura, T., Kodaka, T., Kondo, S., *et al.* Is the cannabinoid CB1 receptor a 2-arachidonoylglycerol receptor? Structural requirements for triggering a Ca²⁺ transient in NG108-15 cells. *J. Biochem.* **122**(4), 890-895 (1997).
2. Kondo, S., Kondo, H., Nakane, S., *et al.* 2-Arachidonoylglycerol, an endogenous cannabinoid receptor agonist: Identification as one of the major species of monoacylglycerols in various rat tissues, and evidence for its generation through Ca²⁺-dependent and -independent mechanisms. *FEBS Lett.* **429**(2), 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, 05/15/2023

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