

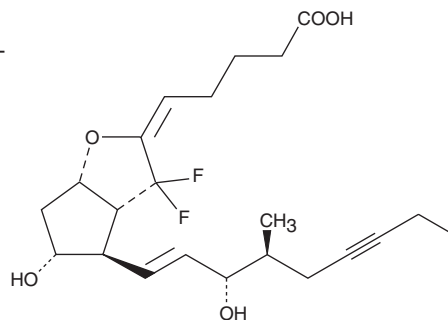
PRODUCT INFORMATION



AFP 07 (free acid)

Item No. 13626

CAS Registry No.: 788799-13-3
Formal Name: 5Z-[(3aR,4R,5R,6aS)-3,3-difluorohexahydro-5-hydroxy-4-[(1E,3S,4S)-3-hydroxy-4-methyl-1-nonen-6-ynyl]-2H-cyclopenta[b]furan-2-ylidene]-pentanoic acid
MF: C₂₂H₃₀F₂O₅
FW: 412.5
Purity: ≥98%
Supplied as: A 5 mg/ml solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

AFP 07 (free acid) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of AFP 07 (free acid) in ethanol and DMF is approximately 30 mg/ml and approximately 25 mg/ml in DMSO.

AFP 07 (free acid) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of AFP 07 (free acid) should be diluted with the aqueous buffer of choice. The solubility of AFP 07 (free acid) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Prostaglandin I₂ is an unstable prostanoid which, through the 'I prostanoid' (IP) receptor, inhibits platelet aggregation and promotes vasodilatation in pulmonary vascular beds. AFP 07 is a 7,7-difluoroprostacyclin derivative that acts as a selective and highly potent agonist for the IP receptor (K_i = 0.561 nM).¹ AFP 07 shows weaker affinity for EP receptors, with K_i values of > 100 nM for EP₁₋₃ and > 10 nM for EP₄.¹ In piglet and rabbit saphenous veins, which express both IP and EP₄, the EP₄ antagonist AH 23848 reduces AFP 07-mediated relaxation, suggesting that AFP 07 activates EP₄ as well as IP, particularly when both receptors are present.²

References

1. Chang, C.-S., Negishi, M., Nakano, T., *et al.* 7,7-Difluoroprostacyclin derivative, AFP-07, a highly selective and potent agonist for the prostacyclin receptor. *Prostaglandins* **53**(2), 83-90 (1997).
2. Jones, R.L. and Chan, K. Distinction between relaxations induced *via* prostanoid EP₄ and IP₁ receptors in pig and rabbit blood vessels. *Br. J. Pharmacol.* **134**(2), 313-324 (2001).

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, 12/08/2025

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