

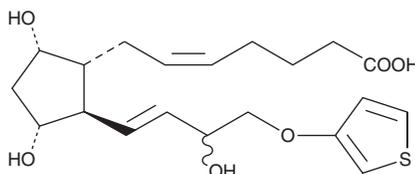
PRODUCT INFORMATION



Tiaprost

Item No. 13632

CAS Registry No.: 71116-82-0
Formal Name: 7-[3,5-dihydroxy-2-[3-hydroxy-4-(3-thienyloxy)-1-buten-1-yl]cyclopentyl]-5-heptenoic acid
Synonym: Iliren
MF: C₂₀H₃₀O₆S
FW: 398.5
Purity: ≥95%
UV/Vis.: λ_{max}: 252 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

Tiaprost 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 tiaprost in these solvents is approximately 30 mg/ml.

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

Description

Prostaglandin F_{2α} plays a central role in luteolysis. Tiaprost is a synthetic analog of PGF_{2α} that is used in veterinary medicine as an estrus-synchronizing agent.¹ It has also been used to treat endometritis and to induce parturition.^{2,3}

References

1. Yadav, V.K., Sudhagar, R.R., and Medhamurthy, R. Apoptosis during spontaneous and prostaglandin F_{2α}-induced luteal regression in the buffalo cow (*Bubalus bubalis*): Involvement of mitogen-activated protein kinases. *Biol. Reprod.* **67**(3), 752-759 (2002).
2. Sheldon, I.M., Lewis, G.S., LeBlanc, S., et al. Defining postpartum uterine disease in cattle. *Theriogenology* **65**(8), 1516-1530 (2006).
3. Holtz, W., Schmidt-Baulain, R., Meyer, H., et al. Control of prostaglandin-induced parturition in sows by injection of the beta-adrenergic blocking agent carazolol or carazolol and oxytocin. *J. Anim. Sci.* **68**(12), 3967-3971 (1990).

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, 05/31/2023

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