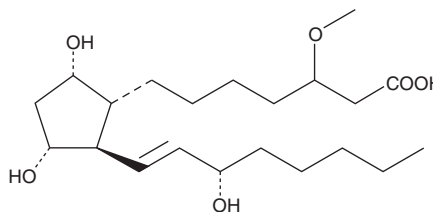


PRODUCT INFORMATION



3-methoxy Prostaglandin F_{1α} Item No. 15170

CAS Registry No.: 54432-43-8
Formal Name: 3-methoxy-9α,11α,15S-trihydroxy-
prost-13Z-en-1-oic acid
Synonym: 3-methoxy PGF_{1α}
MF: C₂₁H₃₈O₆
FW: 386.5
Purity: ≥95%
Supplied as: A 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

3-methoxy Prostaglandin F_{1α} (3-methoxy PGF_{1α}) 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 purged with an inert gas can be used. The solubility of 3-methoxy PGF_{1α} in these solvents is approximately 100, 33, and 23 mg/ml, respectively.

3-methoxy PGF_{1α} is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of 3-methoxy PGF_{1α} should be diluted with the aqueous buffer of choice. The solubility of 3-methoxy PGF_{1α} in PBS (pH 7.2) is approximately 16.7 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

3-methoxy PGF_{1α} is a novel analog of PGF_{1α} with a methoxy group in the 3-position of the upper side chain. There are no published reports describing the pharmacology of 3-methoxy PGF_{1α} at this time.

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, 03/14/2024

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