

PRODUCT INFORMATION



Misoprostol (free acid)-d₅

Item No. 10010333

CAS Registry No.: 1337917-44-8

Formal Name: 9-oxo-11a,16-dihydroxy-16-(methyl-d₃)-prost-13E-en-1-oic-17,17-d₂ acid

MF: C₂₁H₃₁D₅O₅

FW: 373.5

Chemical Purity: ≥98% Misoprostol (free acid)

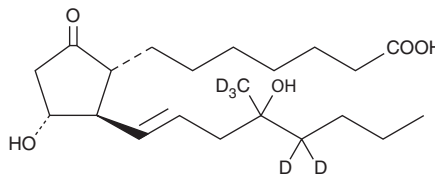
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀

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

Misoprostol (free acid)-d₅ is intended for use as an internal standard for the quantification of misoprostol (free acid) (Item No. 13821) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Misoprostol (free acid)-d₅ 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 misoprostol (free acid)-d₅ in ethanol and DMSO is approximately 50 mg/ml and approximately 100 mg/ml in DMF.

Description

Misoprostol is a PGE₁ analog with agonist activity mediated by EP₂, EP₃, and EP₄ receptors.¹⁻⁴ It has been shown to inhibit the formation of gastric lesions in rats (ED₅₀ = 0.31 μg/kg),² inhibit superoxide generation in human neutrophils (EC₅₀ = 0.35 μM),² and relax fetal rabbit ductus arteriosus (EC₅₀ = 0.36 nM)⁴ in a concentration dependent manner. Misoprostol is commonly used in clinical medicine for the prevention of peptic ulcer disease. It has also been used in conjunction with RU-486 for the oral induction of first trimester abortion. Misoprostol contains a C-1 methyl ester and is readily absorbed and rapidly hydrolyzed in humans to the active free acid.¹

References

1. Walt, R.P. Misoprostol for the treatment of peptic ulcer and anti-inflammatory-drug-induced gastroduodenal ulceration. *N. Engl. J. Med.* **327**, 1575-1580 (1992).
2. Bunce, K.T., Clayton, N.M., Coleman, R.A., et al. GR63799X - a novel prostanoid with selectivity for EP 3 receptors. *Adv. Prostaglandin Thromboxane Leukot. Res.* **21**, 379-382 (1990).
3. Talpain, E., Armstrong, R.A., Coleman, R.A., et al. Characterization of the PGE receptor subtype mediating inhibition of superoxide production in human neutrophils. *Br. J. Pharmacol.* **114**, 1459-1465 (1995).
4. Smith, G.C.S., Coleman, R.A., and McGrath, J.C. Characterization of dilator prostanoid receptors in the fetal rabbit ductus arteriosus. *J. Pharmacol. Exp. Ther.* **271**, 390-396 (1994).

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, 02/26/2024

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