

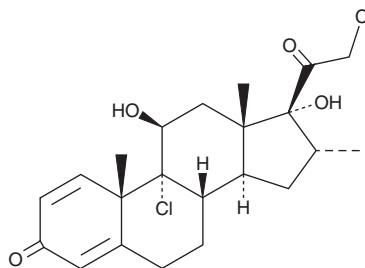
PRODUCT INFORMATION



Mometasone

Item No. 31702

CAS Registry No.: 105102-22-5
Formal Name: (11 β ,16 α)-9,21-dichloro-11,17-dihydroxy-16-methyl-pregna-1,4-diene-3,20-dione
Synonym: (+)-Mometasone
MF: C₂₂H₂₈Cl₂O₄
FW: 427.4
Purity: \geq 90%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Mometasone is supplied as a solid. A stock solution may be made by dissolving the mometasone in the solvent of choice, which should be purged with an inert gas. Mometasone is slightly soluble in methanol.

Description

Mometasone is a synthetic glucocorticoid with anti-inflammatory and anti-allergic activities.¹⁻³ *In vivo*, mometasone (3 mg/kg) reduces pulmonary edema and neutrophil infiltration in a mouse model of chlorine-induced chemical lung injury.¹ Intranasal administration of mometasone (0.2 μ g/animal) reduces nose-rubbing behavior, sneezing, and nasal mucosal levels of IL-4 in a mouse model of allergic rhinitis.² Mometasone also reduces nasal mucosal epithelial destruction, fibrosis, and cilia loss in a rat model of rhinitis medicamentosa induced by oxymetazoline (Item No. 23826).

References

1. Chen, J., Mo, Y., Schlueter, C.F., *et al.* Inhibition of chlorine-induced pulmonary inflammation and edema by mometasone and budesonide. *Toxicol. Appl. Pharmacol.* **272(2)**, 408-413 (2013).
2. Kim, D.H., Kim, B.-Y., Shin, J.-H., *et al.* Intranasal azelastine and mometasone exhibit a synergistic effect on a murine model of allergic rhinitis. *Am. J. Otolaryngol.* **38(2)**, 198-203 (2017).
3. Cam, B., Sari, M., Midi, A., *et al.* Xylitol treats nasal mucosa in rhinitis medicamentosa: An experimental rat model study. *Eur. Arch. Otorhinolaryngol.* **6(11)**, 3123-3130 (2019).

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, 12/08/2022

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