

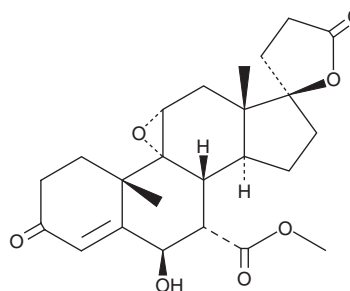
PRODUCT INFORMATION



6 β -hydroxy Eplerenone

Item No. 28859

CAS Registry No.: 209253-80-5
Formal Name: (6 β ,7 α ,11 α ,17 α)-9,11-epoxy-6,17-dihydroxy-3-oxo-pregn-4-ene-7,21-dicarboxylic acid, γ -lactone, methyl ester
MF: C₂₄H₃₀O₇
FW: 430.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

6 β -hydroxy Eplerenone is supplied as a solid. A stock solution may be made by dissolving the 6 β -hydroxy eplerenone in the solvent of choice, which should be purged with an inert gas. 6 β -hydroxy Eplerenone is soluble in methanol.

Description

6 β -hydroxy Eplerenone is a major metabolite of the mineralocorticoid receptor antagonist eplerenone (Item No. 15616).¹ It is formed from eplerenone by the cytochrome P450 (CYP) isoform CYP3A4.

Reference

1. Cook, C.S., Berry, L.M., Kim, D.H., et al. Involvement of CYP3A in the metabolism of eplerenone in humans and dogs: Differential metabolism by CYP3A4 and CYP3A5. *Drug Metab. Dispos.* **30**(12), 1344-1351 (2002).

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