

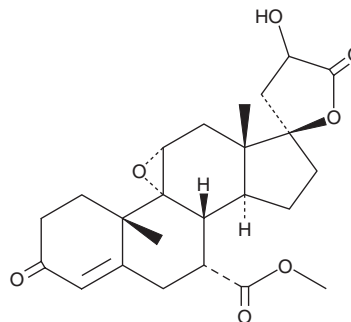
# PRODUCT INFORMATION



## 21-hydroxy Eplerenone

Item No. 28021

**CAS Registry No.:** 334678-67-0  
**Formal Name:** 9,11 $\alpha$ -epoxy-17 $\alpha$ ,21-dihydroxy-3-oxo-pregn-4-ene-7 $\alpha$ ,21-dicarboxylic acid,  $\gamma$ -lactone, methyl ester  
**MF:** C<sub>24</sub>H<sub>30</sub>O<sub>7</sub>  
**FW:** 430.5  
**Purity:**  $\geq$ 98%  
**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

21-hydroxy Eplerenone is supplied as a solid. A stock solution may be made by dissolving the 21-hydroxy Eplerenone in the solvent of choice, which should be purged with an inert gas. 21-hydroxy Eplerenone is slightly soluble in chloroform and methanol.

### Description

21-hydroxy Eplerenone is a major metabolite of the mineralocorticoid receptor antagonist eplerenone (Item No. 15616).<sup>1</sup> 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, 12/19/2022

#### 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