

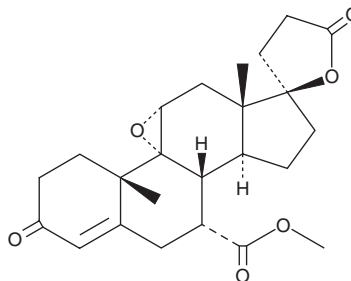
# PRODUCT INFORMATION



## Eplerenone

Item No. 26302

**CAS Registry No.:** 107724-20-9  
**Formal Name:** (7 $\alpha$ ,11 $\alpha$ ,17 $\alpha$ )-9,11-epoxy-17-hydroxy-3-oxo-pregn-4-ene-7,21-dicarboxylic acid,  $\gamma$ -lactone, 7-methyl ester  
**Synonyms:** CGP 30083, (+)-Eplerenone, Epoxymexrenone, SC-6110, SC-66110  
**MF:** C<sub>24</sub>H<sub>30</sub>O<sub>6</sub>  
**FW:** 414.5  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 241 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 5 years



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

### Description

Eplerenone (Item No. 26302) is an analytical reference standard categorized as a diuretic.<sup>1,2</sup> It has been detected as an adverse analytical finding (AAF) during anti-doping testing.<sup>3</sup> This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 15616).

### References

1. Cadwallader, A.B., de la Torre, X., Tieri, A., *et al.* The abuse of diuretics as performance-enhancing drugs and masking agents in sport doping: Pharmacology, toxicology and analysis. *Br. J. Pharmacol.* **161**(1), 1-16 (2010).
2. Reyes, A.J., Leary, W.P., Crippa, G., *et al.* The aldosterone antagonist and facultative diuretic eplerenone: A critical review. *Eur. J. Intern. Med.* **16**(1), 3-11 (2005).
3. World Anti-Doping Agency. *2014 Anti-doping testing figures report* (2014).

#### 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, 07/17/2023

#### 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