

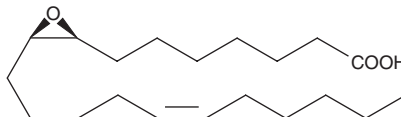
# PRODUCT INFORMATION



## (±)8(9)-EE-14(Z)-E

Item No. 18112

**Formal Name:** 7-((2S,3R)-3-((Z)-undec-5-en-1-yl)oxiran-2-yl)heptanoic acid  
**Synonym:** (±)8,9-Epoxyeicosa-14(Z)-enoic Acid  
**MF:** C<sub>20</sub>H<sub>36</sub>O<sub>3</sub>  
**FW:** 324.5  
**Purity:** ≥95%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



NOTE: Relative stereochemistry shown in chemical structure

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

### Laboratory Procedures

(±)8(9)-EE-14(Z)-E is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of (±)8(9)-EE-14(Z)-E in these solvents is approximately 10 mg/ml.

(±)8(9)-EE-14(Z)-E is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (±)8(9)-EE-14(Z)-E should be diluted with the aqueous buffer of choice. The solubility of (±)8(9)-EE-14(Z)-E in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

(±)8(9)-EE-14(Z)-E is an oxylipin and a metabolite of arachidonic acid (Item Nos. 90010.1 | 90010 | 10006607).<sup>1</sup> (±)8(9)-EE-14(Z)-E increases glomerular membrane albumin permeability and reduces the membrane protective effects of 8(9)-EET in isolated rat glomeruli when used at a concentration of 300 nM.

### Reference

- Sharma, M., McCarthy, E.T., Reddy, D.S., *et al.* 8,9-Epoxyeicosatrienoic acid protects the glomerular filtration barrier. *Prostaglandins Other Lipid Mediat.* **89(1-2)**, 43-51 (2009).

#### 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, 02/02/2024

#### 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