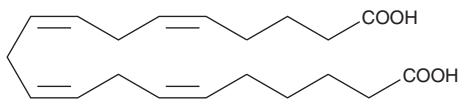


# PRODUCT INFORMATION

## 20-carboxy Arachidonic Acid

Item No. 10007912

**CAS Registry No.:** 79551-84-1  
**Formal Name:** 5Z,8Z,11Z,14Z-eicosatetraenedioic acid  
**Synonyms:** 20-carboxy AA, 20-COOH-AA, FA 20:4;O2  
**MF:** C<sub>20</sub>H<sub>30</sub>O<sub>4</sub>  
**FW:** 334.5  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

20-carboxy Arachidonic acid (20-COOH-AA) 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 20-COOH-AA in these solvents is approximately 100 mg/ml.

20-COOH-AA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 20-COOH-AA should be diluted with the aqueous buffer of choice. The solubility of 20-COOH-AA in 0.1 M Na<sub>2</sub>CO<sub>3</sub> is approximately 1.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

20-COOH-AA is the major metabolite of 20-HETE that is produced in renal tubular epithelial, endothelial, and microvascular smooth muscle cell cultures. This ω-oxidation conversion can take place using purified alcohol dehydrogenases three and four or by microsomes containing recombinant human CYP4F3B.<sup>1</sup> Like 20-HETE, 20-COOH-AA inhibits ion transport in the kidneys. It also produces vasorelaxation of porcine coronary microvessels constricted with endothelin. 20-COOH-AA binds to isolated ligand binding domains of peroxisome proliferator-activated receptor α (PPARα) (K<sub>d</sub> = 0.87 ± 0.12 μM) and PPARγ (K<sub>d</sub> = 1.7 ± 0.5 μM), and is a dual activator of PPARα and PPARγ in a transiently transfected COS-7 cell reporter system.<sup>1</sup>

### Reference

1. Fang, X., Dillon, J.S., Hu, S., *et al.* 20-carboxy-arachidonic acid is a dual activator of peroxisome proliferator-activated receptors α and γ. *Prostaglandins Other Lipid Mediat.* **82(1-4)**, 175-184 (2007).

#### 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, 03/13/2024

#### 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