

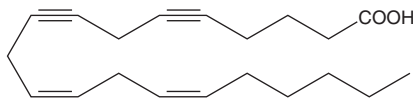
# PRODUCT INFORMATION



## 5,8,11,14-Eicosatetraynoic Acid

Item No. 90120

CAS Registry No.: 1191-85-1  
Synonym: ETYA, FA 20:8  
MF:  $C_{20}H_{24}O_2$   
FW: 296.4  
Purity:  $\geq 98\%$   
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

5,8,11,14-Eicosatetraynoic acid (ETYA) is supplied as a crystalline solid. A stock solution may be made by dissolving the ETYA in the solvent of choice, which should be purged with an inert gas. ETYA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ETYA in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

ETYA is a nonspecific inhibitor of cyclooxygenases (COX) and lipoxygenases (LO).<sup>1</sup> ETYA inhibits human platelet 12-LO and COX-1 with  $IC_{50}$  values of 4  $\mu M$  and 8  $\mu M$ , respectively.<sup>1</sup> ETYA is a PPAR ligand which activates PPAR $\alpha$  and PPAR $\gamma$  chimeras at concentrations of 10  $\mu M$ .<sup>2</sup>

### References

1. Hammarström, S. Selective inhibition of platelet *n*-8 lipoxygenase by 5,8,11-eicosatriynoic acid. *Biochim. Biophys. Acta* **487**, 517-519 (1977).
2. Kliewer, S.A., Lenhard, J.M., Willson, T.M., *et al.* A prostaglandin  $J_2$  metabolite binds peroxisome proliferator-activated receptor  $\gamma$  and promotes adipocyte differentiation. *Cell* **83**, 813-819 (1995).

#### 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, 05/02/2024

#### 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