

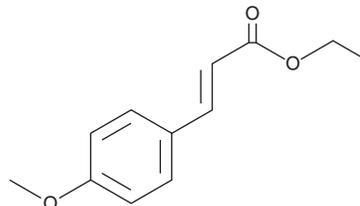
# PRODUCT INFORMATION



## Ethyl *p*-methoxycinnamate

Item No. 11740

**CAS Registry No.:** 24393-56-4  
**Formal Name:** 3-(4-methoxyphenyl)-2E-propenoic acid, ethyl ester  
**Synonyms:** EPMC, Ethyl 4-methoxycinnamate, Ethyl *para*-methoxycinnamate  
**MF:** C<sub>12</sub>H<sub>14</sub>O<sub>3</sub>  
**FW:** 206.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 277, 310 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Hedychium spicatum*



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

### Laboratory Procedures

Ethyl *p*-methoxycinnamate (EPMC) is supplied as a crystalline solid. A stock solution may be made by dissolving the EPMC in the solvent of choice, which should be purged with an inert gas. EPMC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of EPMC in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

EPMC is a cinnamic acid ester that has been found in *K. galanga* and has diverse biological activities.<sup>1-3</sup> It is active against *T. rubrum*, *A. niger*, *S. cerevisiae*, and *E. floccosum* when used at concentrations less than 10 µg/ml.<sup>1</sup> EPMC inhibits COX-1 and COX-2 *in vitro* (IC<sub>50</sub>s = 1.12 and 0.83 µM, respectively).<sup>2</sup> It inhibits microvessel sprouting in isolated rat aortic rings (IC<sub>50</sub> = 91.9 µg/ml).<sup>3</sup> *In vivo*, EPMC (60 mg/kg) reduces IL-1 and TNF-α production and inhibits granuloma formation in a rat model of cotton pellet-induced granuloma formation. It also increases the latency to tail withdrawal in a hot plate test in rats when administered at doses ranging from 200 to 800 mg/kg.

### References

1. Gupta, S.K., Banerjee, A.B., and Achari, B. Isolation of ethyl *p*-methoxycinnamate, the major antifungal principle of *Curcumba zedoaria*. *Lloydia* **39**(4), 218-222 (1976).
2. Umar, M.I., Asmawi, M.Z., Sadikun, A., et al. Bioactivity-guided isolation of ethyl-*p*-methoxycinnamate, an anti-inflammatory constituent, from *Kaempferia galanga* L. extracts. *Molecules* **17**(7), 8720-8734 (2012).
3. Umar, M.I., Asmawi, M.Z., Sadikun, A., et al. Ethyl-*p*-methoxycinnamate isolated from *Kaempferia galanga* inhibits inflammation by suppressing interleukin-1, tumor necrosis factor-α, and angiogenesis by blocking endothelial functions. *Clinics (Sao Paulo)* **69**(2), 134-144 (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, 08/25/2023

#### 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