

# PRODUCT DATA SHEET

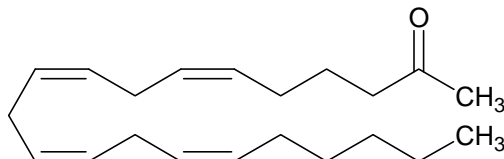


PRODUCT: AACOCH<sub>3</sub>  
Arachidonylmethyl ketone

CATALOG NO.: ST-337

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Colorless oil

MOLECULAR FORMULA: C<sub>21</sub>H<sub>34</sub>O

MOLECULAR WEIGHT: 302.5

PURITY: 98%(TLC)

SOLUBILITY: Soluble in DMSO and ethanol

STORAGE: Store, as supplied, at -20°C for up to 1 year. Store solutions under at atmosphere of argon or nitrogen at -20°C for up to 1 month.

APPLICATION NOTES: Subject to oxidation. Protect from air. Negative control for AACOCF<sub>3</sub><sup>1,2</sup> but stimulates arachidonic acid release *via* possible inhibition of metabolizing or uptake enzymes or *via* direct stimulation of PLA<sub>2</sub>, which has been described<sup>1</sup>.

REFERENCES: 1. A.Street *et al. Biochemistry* 1993 **32** 5935  
2. D.Riendeau *et al. J.Biol.Chem.* 1994 **fo** 15619

**The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.**