

PRODUCT DATA SHEET



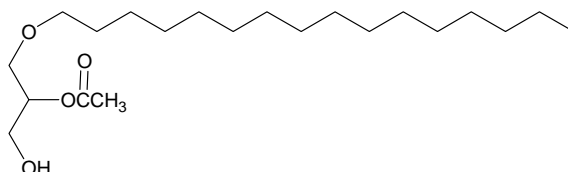
PRODUCT: 1-O-Hexadecyl-2-O-acetyl-sn-glycerol

CAS NO: 77133-35-8

CATALOG NO.: EI-150

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: Waxy solid

MOLECULAR FORMULA: $C_{21}H_{42}O_4$

MOLECULAR WEIGHT: 358.6

PURITY: 95% (TLC) Remainder 1,3-isomer

SOLUBILITY: Soluble in ethanol and acetone

STORAGE: Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 6 months.

APPLICATION NOTES: Avoid acidic or basic conditions. Inhibits protein kinase C activation by diacylglycerols.¹ Stimulates HL-60 cell differentiation.² Biosynthetic precursor of PAF via remodeling and retroconversion/*de novo* routes.³ Induces delayed aggregation of rabbit platelets.⁴ Mitogenic for smooth muscle cells.⁵

- REFERENCES:
1. L.W.Daniel *et al.* *Biochem.Biophys.Res.Comm.* 1988 **151** 291
 2. M.J.C.McNamara *et al.* *Biochem.Biophys.Res.Comm.* 1984 **122** 824
 3. T.-C.Lee *et al.* *J.Biol.Chem.* 1990 **265** 9181
 4. K.Satouchi *et al.* *Arch.Biochem.Biophys.* 1984 **234** 318
 5. L.L.Stoll *et al.* *Cell Regulation* 1989 **1** 13

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.