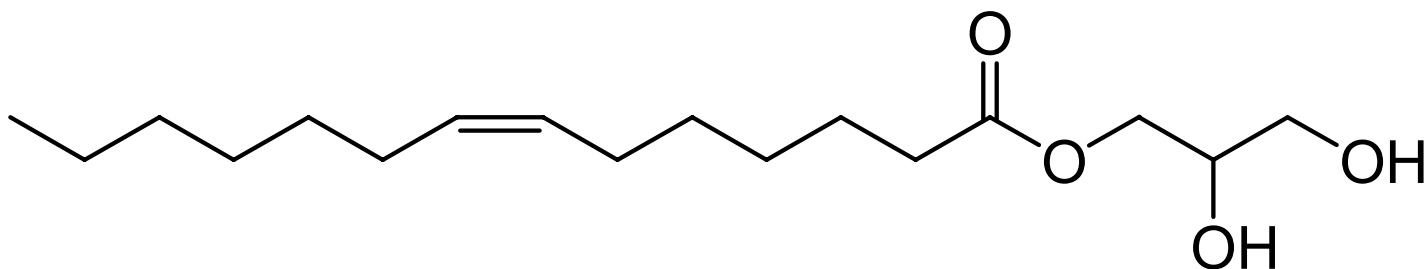


TECHNICAL DATA SHEET

1-(7Z-tetradecenoyl)-rac-glycerol

| | | | |
|---------------------|--|------------------|-----------------|
| Catalog Number | 850530 | Physical state | Oil |
| Purity | 99% | Transition temp. | No data |
| CAS | | CMC | No data |
| Synonyms | 7.7 MAG | pK _a | No data |
| Molec. Formula | C ₁₇ H ₃₂ O ₄ | TLC mobile phase | C:M*, 95:5, v/v |
| MW | 300.434 | Exact Mass | 300.230 |
| Percent composition | C 67.96% H 10.74% O 21.30% | | |
| Stability | Store in <-20°C freezer for up to 6 months | | |
| Solubility | Soluble in ethanol, methanol, water and chloroform | | |
| Web link | 850530 | | |

*C, chloroform; M, methanol



Description:

Crystallizing membrane proteins has been challenging. Recently, lipidic mesophases have been used to isolate the structures of many prokaryotic and eukaryotic membrane proteins, including G-protein receptors (Caffrey and Cherezov, 2009; Caffrey and Porter, 2010). The lipid choice for *in meso* crystallization must be considered (Li, Lee and Caffrey, 2011) because lipid bilayers have different compositions. This methodology has been automated for high-throughput robots (Caffrey and Cherezov, 2009). Avanti Polar Lipids currently offers four modified MAGs to use in *in meso* crystallization studies.

References:

- Caffrey M, Cherezov V (2009) Crystallizing membrane proteins using lipidic mesophases. Nat Protoc. 4(5):706-31
- Caffrey M, Porter C (2010) Crystallizing membrane proteins for structure determination using lipidic mesophases. J Vis Exp.45:1712
- Li D, Lee J, Caffrey M (2011) Crystallizing membrane proteins in lipidic mesophases. A host lipid screen. Cryst Growth Des. 11(2): 530-7

Related products: [Detergents](#)

MSDS: Available at www.avantilipids.com for Product Number 850530