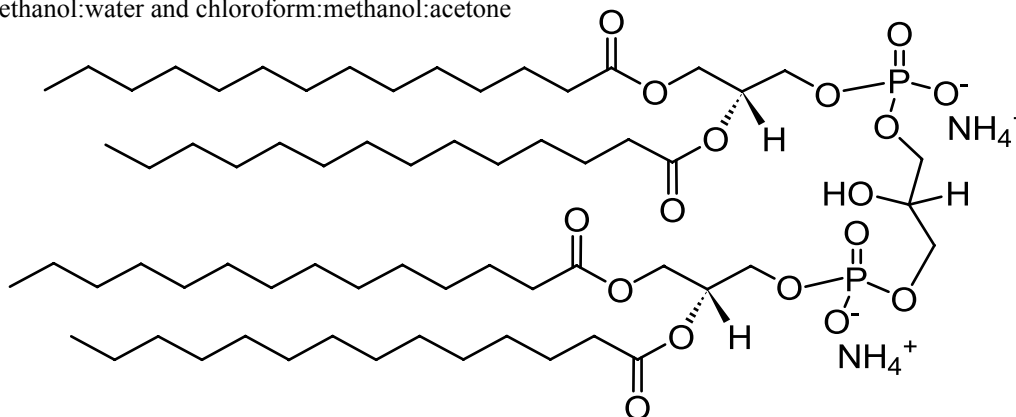


TECHNICAL DATA SHEET

1',3'-bis[1,2-dimyristoyl-*sn*-glycero-3-phospho]-*sn*-glycerol (ammonium salt)

Catalog Number	710332	Physical state	Powder; chloroform solution
Purity	> 99%	Transition temp.	No data
CAS	63988-20-5	CMC	No data
Synonyms	1,1',2,2' Tetramyristoyl Cardiolipin; CL(1'-[14:0/14:0],3'-[14:0/14:0]); 14:0 CA	PKA	No data
Molec. Formula	C ₆₅ H ₁₃₂ N ₂ O ₁₇ P ₂	TLC mobile phase	C:M:W & C:M:A*, 65:25:4, v/v; C:Hexane:M:Acetic Acid 50:30:10:5, v/v, if want to see PG
MW	1,275.690	Exact Mass	1,274.900
Percent composition	C 61.20% H 10.43% N 2.20% O 21.32% P 4.86%		
Stability	Store in <-20°C freezer for 12 months		
Solubility	Soluble in chloroform and alcohol (warmed with water content 3-5%) at concentrations up to 10 mg/mL. Insoluble in water and acetone.		
Web link	710332		

* chloroform:methanol:water and chloroform:methanol:acetone



Description:

Cardiolipin, the mitochondrial-specific phospholipid, helps maintain mitochondrial function, membrane potential and structural support for the inner mitochondrial membrane and the proteins in it (Chicco and Sparagna, 2007). Concentration and composition changes of cardiolipin have been implicated in pathological conditions including ischemia, hypothyroidism, aging, heart failure and cardioskeletal myopathy (Barth syndrome) (Chicco and Sparagna, 2007). Cardiolipin has also been shown as a signaling molecule in apoptosis (Kagan et al, 2006; Ritov et al, 2006).

Product use:

Soluble in chloroform and alcohol (warmed with water content 3-5%) at concentrations up to 10 mg/mL. Insoluble in water and acetone. Also used in liposomes. Please use the following web link for [liposome preparation](#).

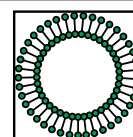
References:

- Huang and Frohman (2009). Lipid Signaling on the Mitochondrial Surface. *Biochim Biophys Acta*. 2009 Jun 17. [Epub ahead of print].
- Chicco and Sparagna (2007). Role of cardiolipin alterations in mitochondrial dysfunction and disease. *Am J Physiol, Cell Physiol* 292(1): C33-44
- Matsumo et al (2006). Lipid domains in bacterial membranes. *Mol Microbiol* 61(5): 1110-7
- Kagan et al (2006). The "pro-apoptotic genes" get out of mitochondria: oxidative lipidomics and redox activity of cytochrome c/cardiolipin complexes. *Chem Biol Interact* 163(1-2): 15-28
- Ritov et al (2006). Analysis of cardiolipin in human muscle biopsy. *J Chromatogr B* 831:63-71

Related products: [cardiolipins](#)

MSDS: see www.avantilipids.com for product number 710332

Avanti Polar Lipids, Inc., 700 Industrial Park Drive, Alabaster, AL 35007-9105, U.S.A.
 Phone • (800) 227-0651 • (205) 663-2494 • Fax (800) 229-1004 • (205) 663-0756
 Email • orders@avantilipids.com • Technical Questions: technical@avantilipids.com
 • Inquiries: info@avantilipids.com • Analytical: analytical@avantilipids.com



Avanti[®]
POLAR LIPIDS, INC.