

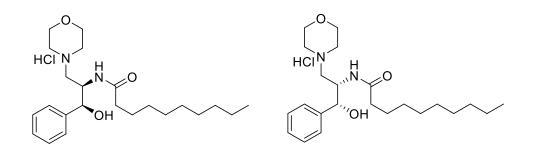
## **PRODUCT DATA SHEET**

## D,L-erythro-PDMP

Catalog number: 1755 Common Name: D,L -*erythro*-1-Phenyl-2decanoylamino-3morpholino-1-propanol • HCl Source: synthetic Solubility: ethanol, methanol, chloroform, DMSO

**CAS number:** 73257-80-4

Molecular Formula: C<sub>23</sub>H<sub>38</sub>N<sub>2</sub>O<sub>3</sub> • HCl Molecular Weight: 427 Storage: -20°C Purity: TLC >98%; HPLC >98%; identity confirmed by MS TLC System: chloroform/methanol/DI water (65:25:4 by vol.) Appearance: solid



## **Application Notes:**

This product inhibits the growth of cells, including cancer cells. D-*threo*-PDMP has been shown to inhibit cell growth by inhibiting the enzyme glucosylceramide synthase<sup>1</sup> but *erythro*-PDMP inhibits growth according to a different mechanism. PDMP has four possible isomers (D-*threo*, L-*threo*, D-*erythro*, and L-*erythro*) due to its two chiral centers. This product (D,L-*erythro*-PDMP) is a mixture of D-*erythro* (1R,2R) and L-*erythro* (1S,2S). The D-*threo* isomer has been shown to be the active glucosyl ceramide synthetase inhibitor.<sup>2</sup> Although *erythro*-PDMP does not inhibit glucosylceramide synthase it does cause cell growth inhibition similar to *threo*-PDMP.<sup>3</sup> This has been suggested as a treatment for cancer.<sup>4</sup> In addition to its stereochemistry, the acyl chain of PDMP has a very marked effect on the intensity of the inhibitory action of the molecule. **Selected References:** 

1. R. Vunnam, N. Radin "Analogs of ceramide that inhibit glucocerebroside synthetase in mouse brain" *Chem Phys Lipids*, Vol. 26(3) pp. 265-278, *1980* 2. N. Radin et al. "Effects of D-*threo*-PDMP, an inhibitor of glucosylceramide synthetase, on expression of cell surface glycolipid antigen and binding to

adhesive proteins by B16 melanoma cells" Journal of Cellular Physiology, Vol. 141(3) pp. 573–583, 1989

3. N. Radin et al. "Effect of an inhibitor of glucosylceramide synthesis on cultured rabbit skin fibroblasts" *Journal of Biochemsitry*, vol. 108:4 pp. 525-530, 1990

4. N. Radin, et al. "Structural and stereochemical studies of potent inhibitors of glucosylceramide synthase and tumor cell growth" *Journal of Lipid Research*, Vol. 36 pp. 611-621, *1995* 

This product is to be used for research only. It is not intended for drug or diagnostic use, human consumption or to be used in food or food additives. Matreya assumes no liability for any use of this product by the end user. We believe the information, offered in good faith, is accurate.