



Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

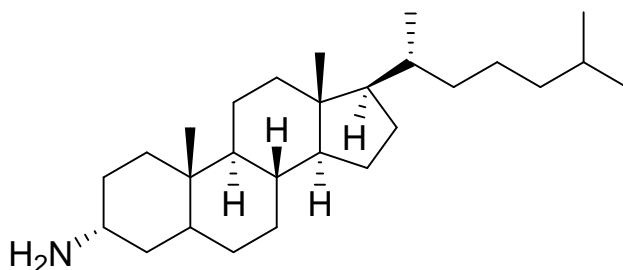
Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name: 3- α -Aminocholestane (3AC)

Product Number: B-0341



Molecular Formula: C₂₇H₄₉N

MW: 387.68

CAS: 2206-20-4

Solubility: EtOH 5 mg/mL

Storage: 4°C. Solutions or suspensions should be stored at -20°C.

Field of Interest: SH2-Domain containing inositol 5-phosphatases (SHIP1 & SHIP2) dephosphorylate the 5-position of PI(3,4,5)P₃ generating PI(3,4)P₂. SHIP2 is ubiquitously expressed while SHIP1 is only found in hematopoietic lineage cells. 3AC is a selective inhibitor of SHIP1 (EC₅₀ = 10 μ M) and shows no inhibition of the other isoform, SHIP2, at concentrations up to 1 mM. 3AC promotes apoptosis of SHIP1-expressing leukemia cells (KG-1) and multiple myeloma cells (OPM) suggesting SHIP1 inhibition is a potential drug target for blood cancers. Mice treated with 3AC show increased numbers of MIR cells in the spleen and lymph nodes and increased numbers of granulocytes.

References: 1) R. Brooks, *et al.*, "SHIP1 Inhibition Increases Immunoregulatory Capacity and Triggers Apoptosis of Hematopoietic Cancer Cells" *J. Immun.*, **2010**, *184*, 3582-3589. 2) G.M. Fuhler, *et al.* "Therapeutic Potential of SH2 Domain-Containing Inositol-5-Phosphatase 1 (SHIP1) and SHIP2 Inhibition in Cancer" *Mol. Med.*, **2012**, *18*, 65-75.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. 3- α -Aminocholestane is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.