



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

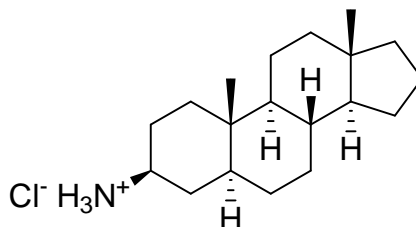
## Technical Data Sheet

### For research use only

Not intended or approved for  
diagnostic or therapeutic use.

**Product Name:** K118 (3 $\beta$ -Amino-5 $\alpha$ -androstanane, hydrochloride salt; SHIP1 Inhibitor)

**Product Number:** B-0344



**Molecular Formula:** C<sub>19</sub>H<sub>34</sub>ClN

**MW:** 311.93

**CAS:** 54156-23-9

**Solubility:** MeOH 5 mg/mL, Water 0.5 mg/mL

**Storage:** 4°C or below

**Field of Interest:** SH2-Domain containing inositol 5-phosphatases (SHIP1 & SHIP2) dephosphorylate the 5-position of PI(3,4,5)P<sub>3</sub> generating PI(3,4)P<sub>2</sub>. SHIP2 is ubiquitously expressed while SHIP1 is found in hematolymphoid cells and mesenchymal stem cells. K118 is a water soluble SHIP1 inhibitor which does not inhibit SHIP2. Treatment with K118 (10 mg/kg) on obese mice on a high fat diet resulted in reduced body mass and lowered fat content when compared to pre-treatment and non-treated control animals. The treated mice had reduced blood glucose and insulin levels in addition to improved glucose tolerance. K118 treatment also reduced the amount of age-related fat accumulation in mice.

**References:** 1) R, Brooks, *et al.* "Coordinate Expansion of Murine Hematopoietic and Mesenchymal Stem Cell Compartments by SHIP1", *Stem Cells*, **2015**, 33, 848-858. doi:10.1002/stem.1902 2) N. Srivastava "A small-molecule inhibitor of SHIP1 reverses age- and diet-associated obesity and metabolic syndrome" *JCI Insight*, **2016**, 1, e88544 doi:10.1172/jci.insight.88544.

**Hazardous Properties and Cautions:** The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. K118 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.