



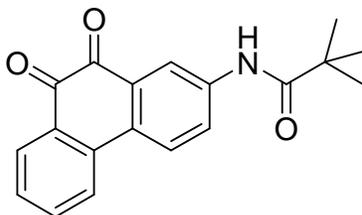
Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Product Name: SF1670 (N-(9,10-Dioxo-9,10-dihydro-phenanthren-2-yl)-2,2-dimethyl-propionamide, PTEN Inhibitor)

Product Number: B-0350



Molecular Formula: C₁₉H₁₇NO₃ **MW:** 307.34 **CAS:** 345630-40-2

Solubility: DMSO (5 mg/mL) or EtOH (1mg/mL, warm)

Storage: 2-8°C

Field of Interest: SF1670 is an inhibitor of PTEN (IC₅₀ = 2 μM). PTEN (Phosphatase and Tensin Homolog deleted on Chromosome 10) is a 3' phosphoinositide phosphatase that converts PI(3,4,5)P₃ to PI(4,5)P₂ thus opposing PKB/Akt activation by PI 3-K. Pretreatment of neutrophils with SF1670 leads to enhanced fMLP-1 induced PI(3,4,5)P₃ signaling effects such as Akt phosphorylation, NADPH oxidase-mediated superoxide formation, cell polarization, chemotactic migration, and phagocytosis. SF1670 also inhibits CD45 (IC₅₀ = 200 nM).

References: 1) Y. Li *et al.* "Pre-treatment with PTEN inhibitor SF1670 augments the efficacy of granulocyte transfusion in a clinically relevant mouse model" *Blood*, **2011**, *117*, 6702-6713.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. SF1670 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.