



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

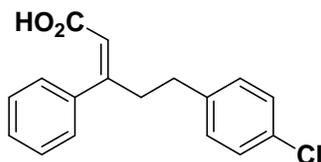
Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name: PS48 (2Z)-5-(4-Chlorophenyl)-3-phenyl-pent-2-enoic acid

Product Number: B-0360



Molecular Formula: C₁₇H₁₅ClO₂

MW: 286.75

CAS: 1180676-32-7

Solubility: DMSO (>20 mg/ml); Ethanol (>10 mg/ml)

Storage: 2~8 °C (protect from light)

Field of Interest: Utilizing low molecular weight compounds to modulate protein kinase activities is a major trend of current pharmaceutical and biological research. PS48 activates phosphoinositide-dependent protein kinase 1 (PDK1) by binding exclusively to PIF-binding pocket rather than the ATP binding site.^[1]

References: 1) A. Stroba *et al.* "3,5-Diphenylpent-2-enoic Acids as Allosteric Activators of the Protein Kinase PDK1: Structure-Activity Relationships and Thermodynamic Characterization of Binding as Paradigms for PIF-Binding Pocket-Targeting Compounds", *J. Med. Chem.*, **2009**, 52 (15), 4683-4693

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. PS48 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.