

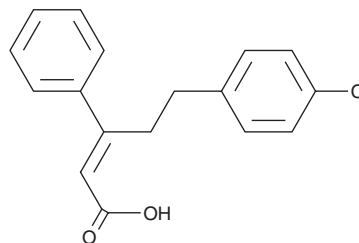
PRODUCT INFORMATION



PS-47

Item No. 41116

CAS Registry No.: 1180676-33-8
Formal Name: 5-(4-chlorophenyl)-3-phenyl-2E-pentenoic acid
MF: C₁₇H₁₅ClO₂
FW: 286.8
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PS-47 is supplied as a solid. A stock solution may be made by dissolving the PS-47 in the solvent of choice, which should be purged with an inert gas. PS-47 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

PS-47 is an inactive derivative of the 3-phosphoinositide-dependent protein kinase 1 (PDK1) allosteric activator PS-48 that has been used as a negative control for the kinase activity of PDK1.^{1,2}

References

1. Stroba, A., Schaeffer, F., Hindie, V., 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.* **52(15)**, 4683-4693 (2009).
2. Querfurth, H., Marshall, J., Parang, K., et al. A PDK-1 allosteric agonist neutralizes insulin signaling derangements and beta-amyloid toxicity in neuronal cells and in vitro. *PLoS One* **17(1)**, e0261696 (2022).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/24/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM