

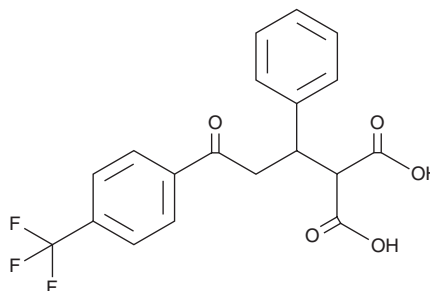
PRODUCT INFORMATION



PS-210

Item No. 30923

CAS Registry No.: 1221962-86-2
Formal Name: 2-[3-oxo-1-phenyl-3-[4-(trifluoromethyl)phenyl]propyl]-propanedioic acid
MF: C₁₉H₁₅F₃O₅
FW: 380.3
Purity: ≥98%
UV/Vis.: λ_{max}: 235 nm
Supplied as: A crystalline 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-210 is supplied as a crystalline solid. A stock solution may be made by dissolving the PS-210 in the solvent of choice, which should be purged with an inert gas. PS-210 is soluble in DMSO.

Description

PS-210 is an activator of phosphoinositide-dependent protein kinase 1 (PDK1) that increases its thermal stability.^{1,2} It is selective for PDK-1 over a panel of 120 kinases at 10 μM.¹ PS-210 (2 μM) increases the activity of wild-type PDK1 but not PDK1 with a mutation in the PIF-binding pocket.

References

1. Busschots, K., Lopez-Garcia, L.A., Lammi, C., *et al.* Substrate-selective inhibition of protein kinase PDK1 by small compounds that bind to the PIF-pocket allosteric docking site. *Chem. Biol.* **19(9)**, 1152-1163 (2012).
2. Wilhem, A., Lopez-Garcia, L.A., Busschots, K., *et al.* 2-(3-Oxo-1,3-diphenylpropyl)malonic acids as potent allosteric ligands of the PIF pocket of phosphoinositide-dependent kinase-1: Development and prodrug concept. *J. Med. Chem.* **55(22)**, 9817-9830 (2012).

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, 12/12/2022

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