

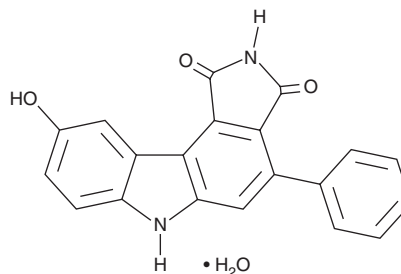
PRODUCT INFORMATION



PD 407824 (hydrate)

Item No. 45243

CAS Registry No.: 1177150-89-8
Formal Name: 9-hydroxy-4-phenyl-pyrrolo[3,4-c]carbazole-1,3(2H,6H)-dione, monohydrate
MF: C₂₀H₁₂N₂O₃ • H₂O
FW: 346.3
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

PD 407824 (hydrate) is supplied as a solid. A stock solution may be made by dissolving the PD 407824 (hydrate) in the solvent of choice, which should be purged with an inert gas. PD 407824 (hydrate) is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

PD 407824 is an inhibitor of the checkpoint kinases Chk1 and WEE1 (IC₅₀s = 47 and 97 nM, respectively).¹ It is selective for Chk1 and WEE1 over PKC (IC₅₀ = 3.4 μM) and Cdk4 (IC₅₀ = 3.75 μM), as well as c-Src and the PDGF and FGF receptors (IC₅₀s = >50 μM for all), and other CDKs (IC₅₀s = >50 μM).² PD 407824 sensitizes SKOV3 and OVCAR-3 ovarian cancer cells, as well as cisplatin-resistant A2780cis cells, to cisplatin when used at a concentration of 0.5 μM.³ It also sensitizes C2C12 myoblasts to bone morphogenetic protein 4 (BMP4) and, when used in combination with BMP4, inhibits myotube formation and induces myoblasts to differentiate into mature osteoblasts.⁴ PD 407824, in combination with BMP4, induces human embryonic stem cells (hESCs) to differentiate into cells with mesoderm or cytotrophoblast stem cell lineages.

References

1. Smaill, J.B., Baker, E.N., Booth, R.J., et al. Synthesis and structure-activity relationships of N-6 substituted analogues of 9-hydroxy-4-phenylpyrrolo[3,4-c]carbazole-1,3(2H,6H)-diones as inhibitors of Wee1 and Chk1 checkpoint kinases. *Eur. J. Med. Chem.* **43(6)**, 1276-1296 (2008).
2. Palmer, B.D., Thompson, A.M., Booth, R.J., et al. 4-Phenylpyrrolo[3,4-c]carbazole-1,3(2H,6H)-dione inhibitors of the checkpoint kinase Wee1. Structure-activity relationships for chromophore modification and phenyl ring substitution. *J. Med. Chem.* **49(16)**, 4896-4911 (2006).
3. Arora, S., Bisanz, K.M., Peralta, L.A., et al. RNAi screening of the kinome identifies modulators of cisplatin response in ovarian cancer cells. *Gynecol. Oncol.* **118(3)**, 220-227 (2010).
4. Feng, L., Cook, B., Tsai, S.-Y., et al. Discovery of a small-molecule BMP sensitizer for human embryonic stem cell differentiation. *Cell Rep.* **15(9)**, 2063-2075 (2016).

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, 02/17/2026

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