

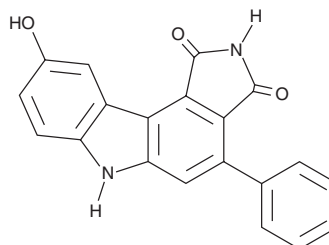
PRODUCT INFORMATION



PD 407824

Item No. 25989

CAS Registry No.: 622864-54-4
Formal Name: 9-hydroxy-4-phenyl-pyrrolo[3,4-c]carbazole-1,3(2H,6H)-dione
MF: C₂₀H₁₂N₂O₃
FW: 328.3
Purity: ≥98%
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 is supplied as a solid. A stock solution may be made by dissolving the PD 407824 in the solvent of choice. PD 407824 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 18 mg/ml.

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), Cdk4 (IC₅₀ = 3.75 μM), as well as c-Src and the PDGF and FGF receptors (IC₅₀s = >50 μM for all), and other Cdk (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 morphogenic 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 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, 03/09/2023

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