

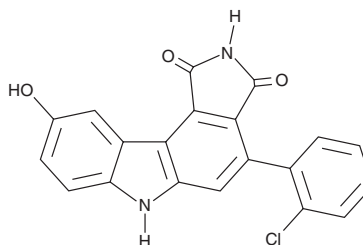
PRODUCT INFORMATION



Wee1 Inhibitor

Item No. 35391

CAS Registry No.: 622855-37-2
Formal Name: 4-(2-chlorophenyl)-9-hydroxy-pyrrolo[3,4-c]carbazole-1,3(2H,6H)-dione
MF: $C_{20}H_{11}ClN_2O_3$
FW: 362.8
Purity: $\geq 95\%$
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

Wee1 inhibitor is supplied as a solid. A stock solution may be made by dissolving the Wee1 inhibitor in the solvent of choice, which should be purged with an inert gas. Wee1 inhibitor is soluble in organic solvents such as ethanol and DMSO. The solubility of Wee1 inhibitor in these solvents is approximately 10 mg/ml.

Description

Wee1 inhibitor is an inhibitor of the checkpoint kinase Wee1 ($IC_{50} = 11$ nM).¹ It is selective for Wee1 over checkpoint kinase 1 (Chk1; $IC_{50} = 440$ nM). Wee1 inhibitor also inhibits the ATPase activity of LptB, an enzyme involved in the transport and assembly of LPS ($IC_{50} = 25$ μM for the *E. coli* enzyme).²

References

1. 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).
2. Gronenberg, L.S. and Kahne, D. Development of an activity assay for discovery of inhibitors of lipopolysaccharide transport. *J. Am. Chem. Soc.* **132**(8), 2518-2519 (2010).

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