

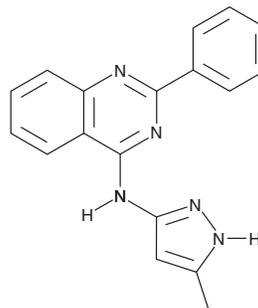
# PRODUCT INFORMATION



## CAY10796

Item No. 21190

**CAS Registry No.:** 404828-14-4  
**Formal Name:** N-(5-methyl-1H-pyrazol-3-yl)-2-phenyl-4-quinazolinamine  
**MF:** C<sub>18</sub>H<sub>15</sub>N<sub>5</sub>  
**FW:** 301.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 210, 249, 331 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

CAY10796 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10796 in the solvent of choice. CAY10796 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of CAY10796 in these solvents is approximately 0.15 and 0.25 mg/ml, respectively.

### Description

CAY10796 is an aminopyrazole ATP-competitive inhibitor of GSK3, with 34% inhibition when used at a concentration of 2.5 μM.<sup>1</sup> It inhibits androgen receptor transactivation in 22Rv1, LNCaP, and LNCaP-SSR cell lines in a dose-dependent manner. It promotes nuclear export of the androgen receptor and decreases translocation to the nucleus in PC3 and PCa prostate cancer cells, respectively.<sup>1,2</sup> In HEK293 cells expressing the rat Na<sub>v</sub>1.2 channel, pretreatment with CAY10796 dose-dependently potentiates peak current densities.<sup>3</sup>

### References

1. Rinnab, L., Schütz, S.V., Diesch, J., *et al.* Inhibition of glycogen synthase kinase-3 in androgen-responsive prostate cancer cell lines: Are GSK inhibitors therapeutically useful? *Neoplasia* **10(6)**, 624-634 (2008).
2. Schütz, S.V., Cronauer, M.V., and Rinnab, L. Inhibition of glycogen synthase kinase-3β promotes nuclear export of the androgen receptor through a CRM1-dependent mechanism in prostate cancer cell lines. *J. Cell. Biochem.* **109(6)**, 1192-1200 (2010).
3. James, T.F., Nenov, M.N., Wildburger, N.C., *et al.* The Na<sub>v</sub>1.2 channel is regulated by GSK3. *Biochim. Biophys. Acta* **1850(4)**, 832-844 (2015).

#### 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, 02/07/2023

#### 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