

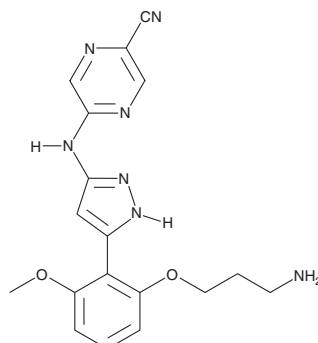
# PRODUCT INFORMATION



**LY2606368**

Item No. 21490

**CAS Registry No.:** 1234015-52-1  
**Formal Name:** 5-[[5-[2-(3-aminopropoxy)-6-methoxyphenyl]-1H-pyrazol-3-yl]amino]-2-pyrazinecarbonitrile  
**Synonym:** Prexasertib  
**MF:** C<sub>18</sub>H<sub>19</sub>N<sub>7</sub>O<sub>2</sub>  
**FW:** 365.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 214, 295 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.

## Description

LY2606368 is a checkpoint kinase 1 (Chk1) inhibitor with a K<sub>i</sub> value of 0.9 nM against the purified target and IC<sub>50</sub> values of <13 nM in a viability study of multiple colorectal cancer cell lines.<sup>1,2</sup> Inhibition of Chk1 causes double-strand DNA breakage in cells leading to an excessive cell damage burden and subsequent cell death.<sup>1,2</sup> *In vitro*, LY2606368 inhibits the doxorubicin-activated G<sub>2</sub>/M checkpoint in p53-deficient HeLa cells with an EC<sub>50</sub> value of 9 nM.<sup>1</sup> LY2606368, at 25 nM, also significantly induces apoptosis and inhibits colony formation in AGS and MKN1 gastric cancer cells.<sup>3</sup>

## References

1. King, C., Diaz, H.B., McNeely, S., *et al.* LY2606368 causes replication catastrophe and antitumor effects through CHK1-dependent mechanisms. *Mol. Cancer Ther.* **14**(9), 2004-2013 (2015).
2. Manic, G., Signore, M., Sistigu, A., *et al.* CHK1-targeted therapy to deplete DNA replication-stressed, p53-deficient, hyperdiploid colorectal cancer stem cells. *Gut* **67**(5), 903-917 (2017).
3. Yin, Y., Shen, Q., Zhang, P., *et al.* Chk1 inhibition potentiates the therapeutic efficacy of PARP inhibitor BMN673 in gastric cancer. *Am. J. Cancer Res.* **7**(3), 473-483 (2017).

### 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, 12/15/2022

## 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