

# PRODUCT INFORMATION

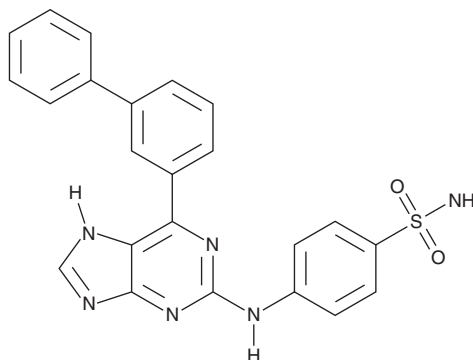


## CDK2 Inhibitor 73

Item No. 42335

**CAS Registry No.:** 2079895-42-2  
**Formal Name:** 4-[[6-[1,1'-biphenyl]-3-yl]-9H-purin-2-yl]amino-benzenesulfonamide  
**Synonyms:** CDK2-IN-4, CDK2-IN-73, CDK2 Inhibitor 4, Cyclin-Dependent Kinase 2-IN-4, Cyclin-Dependent Kinase 2-IN-73, Cyclin-Dependent Kinase 2 Inhibitor 4, Cyclin-Dependent Kinase 2 Inhibitor 73

**MF:** C<sub>23</sub>H<sub>18</sub>N<sub>6</sub>O<sub>2</sub>S  
**FW:** 442.5  
**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

CDK2 inhibitor 73 is supplied as a solid. A stock solution may be made by dissolving the CDK2 inhibitor 73 in the solvent of choice, which should be purged with an inert gas. CDK2 inhibitor 73 is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

CDK2 inhibitor 73 is an inhibitor of cyclin-dependent kinase 2 (Cdk2; IC<sub>50</sub> = 0.044 μM).<sup>1</sup> It is selective for Cdk2 over Cdk1, Cdk4, Cdk7, and Cdk9 at 100 μM.

### Reference

1. Coxon, C.R., Anscombe, E., Harnor, S.J., *et al.* Cyclin-dependent kinase (CDK) inhibitors: Structure-activity relationships and insights into the CDK-2 selectivity of 6-substituted 2-arylaminopurines. *J. Med. Chem.* **60**(5), 1746-1767 (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, 11/14/2024

#### 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