

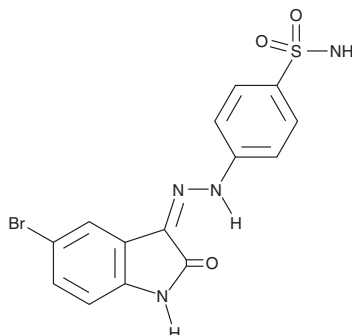
# PRODUCT INFORMATION



## Cdk2 Inhibitor II

Item No. 15154

**CAS Registry No.:** 222035-13-4  
**Formal Name:** 4-[2-(5-bromo-1,2-dihydro-2-oxo-3H-indol-3-ylidene)hydrazinyl]-benzenesulfonamide  
**Synonyms:** Cyclin-dependent kinase 2 Inhibitor II, SC-221409  
**MF:** C<sub>14</sub>H<sub>11</sub>BrN<sub>4</sub>O<sub>3</sub>S  
**FW:** 395.2  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 213, 257, 395 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

Cdk2 inhibitor II is supplied as a crystalline solid. A stock solution may be made by dissolving the Cdk2 inhibitor II in the solvent of choice, which should be purged with an inert gas. Cdk2 inhibitor II is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

### Description

The cyclin-dependent kinase 2 (Cdk2) works with cyclins A or E to regulate S phase and G<sub>2</sub>-M transition during the cell cycle.<sup>1</sup> Cdk2 inhibitor II is a 3-(benzylidene)indolin-2-one analog that selectively and potently inhibits Cdk2 (IC<sub>50</sub> = 60 nM).<sup>2</sup> It is cell permeable, reversible, and ATP-competitive.<sup>2</sup>

### References

1. Funk, J.O. Cell cycle checkpoint genes and cancer. *Encyclopedia of life sciences*. John Wiley & Sons (2005).
2. Davis, S.T., Benson, B.G., Bramson, H.N., *et al.* Prevention of chemotherapy-induced alopecia in rats by CDK inhibitors. *Science* **291(5501)**, 134-137 (2001).

#### 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, 11/14/2022

#### 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