

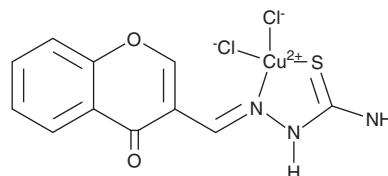
# PRODUCT INFORMATION



## Akt Inhibitor XI

Item No. 18604

**CAS Registry No.:** 902779-59-3  
**Formal Name:** (SP-4-3)-dichloro[(2Z)-2-[(4-oxo-4H-1-benzopyran-3-yl)methylene]hydrazinecarbothioamide-κN<sup>2</sup>,κS]-copper  
**Synonym:** FPA-124  
**MF:** C<sub>11</sub>H<sub>9</sub>Cl<sub>2</sub>CuN<sub>3</sub>O<sub>2</sub>S  
**FW:** 381.7  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 273, 308 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

Akt inhibitor XI is supplied as a crystalline solid. A stock solution may be made by dissolving the Akt inhibitor XI in the solvent of choice, which should be purged with an inert gas. Akt inhibitor XI is soluble in the organic solvent DMSO at a concentration of approximately 0.1 mg/ml.

### Description

The kinase Akt (also known as protein kinase B or PKB) modulates cell proliferation, metabolism, and survival, as well as angiogenesis.<sup>1,2</sup> Akt inhibitor XI is a cell-permeable, copper-containing 3-formylchromone derivative that inhibits Akt in an array of cancer cells (IC<sub>50</sub>s = 10-34 μM).<sup>3</sup> It also causes NF-κB inactivation in an orthotopic pancreatic tumor model using COLO 357 cells.<sup>3</sup> Molecular modeling indicates that this inhibitor interacts with the pleckstrin homology and kinase domains of Akt. Akt inhibitor XI is commonly used in the range of 1-20 μM to assess the role of Akt in cellular responses.<sup>4-6</sup>

### References

1. Manning, B.D. and Cantley, L.C. AKT/PKB signaling: Navigating downstream. *Cell* **129**(7), 1261-1274 (2007).
2. Yuan, T.L. and Cantley, L.C. PI3K pathway alterations in cancer: Variations on a theme. *Oncogene* **27**(41), 5497-5510 (2008).
3. Barve, V., Ahmed, F., Adsule, S., et al. Synthesis, molecular characterization, and biological activity of novel synthetic derivatives of chromen-4-one in human cancer cells. *J. Med. Chem.* **49**(13), 3800-3808 (2006).
4. Frampton, G., Invernizzi, P., Bernuzzi, F., et al. Interleukin-6-driven progranulin expression increases cholangiocarcinoma growth by an Akt-dependent mechanism. *Gut* **61**(2), 268-277 (2012).
5. Rybchyn, M.S., Slater, M., Conlgrave, A.D., et al. An Akt-dependent increase in canonical Wnt signaling and a decrease in sclerostin protein levels are involved in strontium ranelate-induced osteogenic effects in human osteoblasts. *J. Biol. Chem.* **286**(27), 23771-23779 (2011).
6. Zareen, N., Biswas, S.C., and Greene, L.A. A feed-forward loop involving Trib3, Akt and FoxO mediates death of NGF-deprived neurons. *Cell Death Differ.* **20**(12), 1719-1730 (2013).

#### 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, 11/04/2022

#### 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