

# PRODUCT INFORMATION

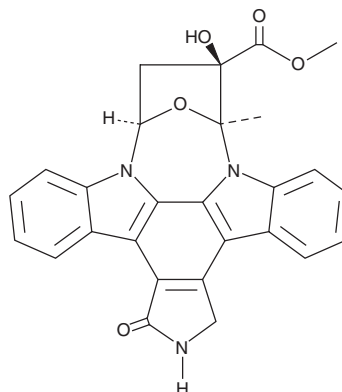


## K252a

Item No. 11338

**CAS Registry No.:** 99533-80-9  
**Formal Name:** 2,3,9S,10R,11,12R-hexahydro-10-hydroxy-9-methyl-1-oxo-9,12-epoxy-1H-diindolo[1,2,3-fg:3',2',1'-kl]pyrrolo[3,4-i][1,6]benzodiazocine-10-carboxylic acid, methyl ester

**Synonym:** SF 2370  
**MF:** C<sub>27</sub>H<sub>21</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 467.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 290 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

K252a is supplied as a crystalline solid. A stock solution may be made by dissolving the K252a in the solvent of choice, which should be purged with an inert gas. K252a is soluble in organic solvents such as methanol, DMSO, and dichloromethane. The solubility of K252a in these solvents is approximately 2, 50, and 5 mg/ml, respectively.

### Description

K252a is a staurosporine analog isolated from *Nocardiosis* sp. soil fungi that inhibits protein kinase (PK) C, PKA, Ca<sup>2+</sup>/calmodulin-dependent kinase type II, and phosphorylase kinase with IC<sub>50</sub> values of 470, 140, 270, and 1.7 nM, respectively.<sup>1,2</sup> Because it inhibits neurotrophin receptor tyrosine kinases, K252a at 100-500 nM has been used to suppress trophoblast proliferation and increase apoptosis associated with the disruption of mitochondrial functions in cultured choriocarcinoma cells.<sup>3</sup> Recently, K252a has been shown to inhibit PRK1 (IC<sub>50</sub> = 3.2 nM *in vitro*), a PKC-related kinase that phosphorylates histone H3 at threonine 11 and is involved in androgen-dependent gene expression.<sup>4</sup>

### References

1. Yasuzawa, T., Iida, T., Yoshida, M., *et al.* The structures of the novel protein kinase C inhibitors K-252a, b, c and d. *J. Antibiot. (Tokyo)* **39**(8), 1072-1078 (1986).
2. Davis, P.D., Hill, C.H., Lawton, G., *et al.* Inhibitors of protein kinase C. 1.1 2,3-bisarylmales. *J. Med. Chem.* **35**(1), 177-184 (1992).
3. Kawamura, N., Kawamura, K., Manabe, M., *et al.* Inhibition of brain-derived neurotrophic factor/tyrosine kinase B signaling suppresses choriocarcinoma cell growth. *Endocrinology* **151**(7), 3006-3014 (2010).
4. Köhler, J., Erlenkamp, G., Eberlin, A., *et al.* Lestaurtinib inhibits histone phosphorylation and androgen-dependent gene expression in prostate cancer cells. *PLoS One* **7**(4), (2012).

#### 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, 08/16/2023

#### 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](http://WWW.CAYMANCHEM.COM)