

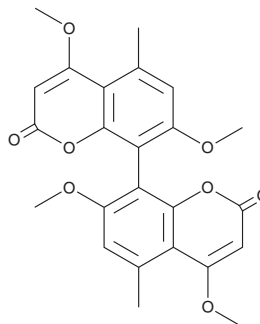
# PRODUCT INFORMATION



## Kotanin A

Item No. 24917

**CAS Registry No.:** 27909-08-6  
**Formal Name:** (8S)- 4,4',7,7'-tetramethoxy-5,5'-dimethyl-[8,8'-bi-2H-1-benzopyran]-2,2'-dione  
**Synonym:** (+)-Kotanin  
**MF:** C<sub>24</sub>H<sub>22</sub>O<sub>8</sub>  
**FW:** 438.4  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungi/*Aspergillus* sp.



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

### Laboratory Procedures

Kotanin A is supplied as a solid. A stock solution may be made by dissolving the kotanin A in the solvent of choice. Kotanin A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

### Description

Kotanin A is a fungal metabolite produced by *A. glaucus* (*A. clavatus*), *A. alliaceus*, and *A. niger*.<sup>1,2</sup> It is toxic to day-old chicks (LD<sub>100</sub> = 62.5 mg/kg) but has no effect in rats.<sup>1</sup>

### References

1. Buchi, G., Klaubert, D.H., Shank, R.C., *et al.* Structure and synthesis of kotanin and desmethylkotanin, metabolites of *Aspergillus glaucus*. *J. Org. Chem.* **36**(8), 1143-1147 (1971).
2. Cutler, H.G., Crumley, F.G., Cox, R.H., *et al.* Orlandin: A nontoxic fungal metabolite with plant growth inhibiting properties. *J. Agric. Food Chem.* **27**(3), 592-595 (1979).

#### 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, 12/19/2022

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