

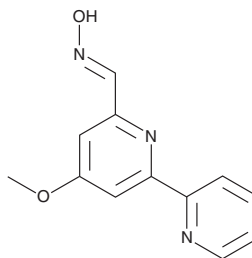
# PRODUCT INFORMATION



## Caerulomycin A

Item No. 27625

**CAS Registry No.:** 21802-37-9  
**Formal Name:** (E)-4-methoxy-[2,2'-bipyridine]-6-carboxaldehyde oxime  
**MF:** C<sub>12</sub>H<sub>11</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 229.2  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Bacterium/*Streptomyces* sp.



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

### Laboratory Procedures

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

Caerulomycin A is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Caerulomycin A is a fungal metabolite originally isolated from *Actinoalloteichus* with antifungal and immunosuppressant activities.<sup>1-4</sup> It inhibits the growth of *C. albicans*, *C. glabrata*, and *C. krusei* (MICs = 0.78-1.56, 0.39-0.78, and 0.78-1.56 µg/ml, respectively).<sup>1</sup> Caerulomycin A induces expansion of CD4<sup>+</sup>Foxp3<sup>+</sup> regulatory T cells (Tregs) and decreases the number of Th1 and T17 cells *in vitro* via increased TGF-β-mediated Smad3 activity and reduces IFN-γ-induced STAT1 signaling.<sup>2</sup> *In vivo*, caerulomycin A (10 mg/kg) reduces IL-6, TNF-α, and IFN-γ production, inflammation, and synovitis in a mouse model of collagen-induced arthritis. Caerulomycin A suppresses the differentiation of Th2 cells and reduces levels of IL-4, IL-5, IL-13, and IgE and eosinophil lung infiltration in a mouse model of ovalbumin-induced asthma.<sup>3</sup> It also increases production of Tregs, reduces production of Th1, Th17, and CD8 T cells, and reduces disease severity in a mouse model of experimental autoimmune encephalomyelitis (EAE).<sup>4</sup>

### References

1. Ambavane, V., Tokdar, P., Parab, R., *et al.* Caerulomycin A - An antifungal compound isolated from marine actinomycetes. *Adv. Microbiol.* **4(9)**, 567-578 (2014).
2. Gurram, R.K., Kujur, W., Maurya, S.K., *et al.* Caerulomycin A enhances transforming growth factor-β (TGF-β)-Smad3 protein signaling by suppressing interferon-γ (IFN-γ)-signal transducer and activator of transcription 1 (STAT1) protein signaling to expand regulatory T cells (Tregs). *J. Biol. Chem.* **289(25)**, 17515-17528 (2014).
3. Kujur, W., Gurram, R.K., Haleem, N., *et al.* Caerulomycin A inhibits Th2 cell activity: A possible role in the management of asthma. *Sci. Rep.* **5**, 15396 (2015).
4. Kujur, W., Gurram, R.K., Maurya, S.K., *et al.* Caerulomycin A suppresses the differentiation of naïve T cells and alleviates the symptoms of experimental autoimmune encephalomyelitis. *Autoimmunity* **50(5)**, 317-328 (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, 12/22/2022

#### 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