

PRODUCT INFORMATION



Acetylaszonalenin

Item No. 25207

CAS Registry No.: 42230-55-7
Formal Name: (5aS,13aR,14aS)-5-acetyl-14a-(1,1-dimethyl-2-propen-1-yl)-5a,13a,14,14a-tetrahydro-indolo[3',2':4,5]pyrrolo[2,1-c][1,4]benzodiazepine-7,13(5H,12H)-dione

MF: C₂₅H₂₅N₃O₃

FW: 415.5

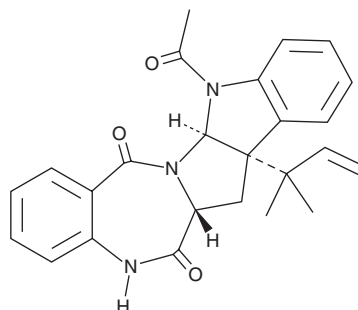
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Neosartorya spinosa*



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

Laboratory Procedures

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

Description

Acetylaszonalenin is a fungal metabolite that has been found in *A. flavipes*.¹ It acts as an antagonist of neurokinin-1 (NK₁) receptors, inhibiting binding of substance P (Item No. 24035) with a K_i value of 170 μM.

Reference

1. Barrow, C.J. and Sun, H.H. Spiroquinazoline, a novel substance P inhibitor with a new carbon skeleton, isolated from *Aspergillus flavipes*. *J. Nat. Prod.* **57(4)**, 471-476 (1994).

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, 08/23/2022

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