

PRODUCT INFORMATION



DSDP

Item No. 30156

CAS Registry No.: 1777807-64-3
Formal Name: (5'aS,10'aS)-1',5'a,6',10'a-tetrahydro-dispiro[2H-indene-2,3'-[3H,5H,8H,10H]bisthiazolo[3,4-a:3',4'-d]pyrazine-8',2''-[2H]indene]-1,1'',3,3'',5',10'-hexone

Synonym: ZINC03129319

MF: C₂₄H₁₄N₂O₆S₂

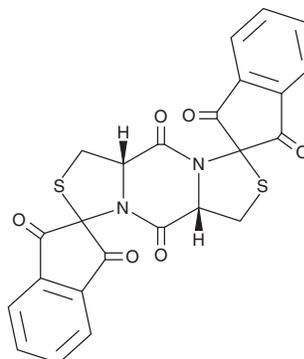
FW: 490.5

Purity: ≥98%

Supplied as: A 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

DSDP is supplied as a solid. A stock solution may be made by dissolving the DSDP in the solvent of choice, which should be purged with an inert gas. DSDP is soluble in the organic solvent DMSO.

Description

DSDP is an activator of stimulator of interferon genes (STING).¹ It induces expression of an IFN-stimulated gene 54 (ISG54) luciferase reporter in HepAD38 cells that constitutively express human cyclic GMP-AMP synthase (cGAS) and STING. DSDP (50 μM) increases *IFNB1*, *TNFA*, and *IL29* expression in HepG2 cells expressing wild-type human, but not C-terminal truncated human or mouse, STING, as well as in isolated human peripheral blood mononuclear cells (PBMCs). It inhibits yellow fever, dengue, and Zika virus replication in THF fibroblasts at the same concentration.

Reference

1. Liu, B., Tang, L., Zhang, X., *et al.* A cell-based high throughput screening assay for the discovery of cGAS-STING pathway agonists. *Antiviral Res.* **147**, 37-46 (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, 11/07/2022

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