

# PRODUCT INFORMATION



**DS-3032**

Item No. 41838

**CAS Registry No.:** 1398568-47-2

**Formal Name:** 2,6-anhydro-5-[[[(3'R,4'S,5'R)-6''-chloro-4'-(2-chloro-3-fluoro-4-pyridinyl)-1'',2''-dihydro-4,4-dimethyl-2''-oxodispiro[cyclohexane-1,2'-pyrrolidine-3',3''-[3H]indol]-5'-yl]carbonyl]amino]-3,4,5-trideoxy-L-erythro-hexonamide

**Synonyms:** DS-3032b, Milademetan, RAIN-32

**MF:** C<sub>30</sub>H<sub>34</sub>Cl<sub>2</sub>FN<sub>5</sub>O<sub>4</sub>

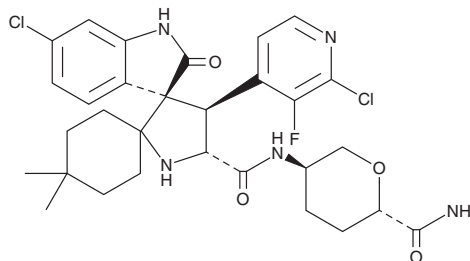
**FW:** 618.5

**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

DS-3032 is supplied as a solid. A stock solution may be made by dissolving the DS-3032 in the solvent of choice, which should be purged with an inert gas. DS-3032 is sparingly soluble (1-10 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

## Description

DS-3032 is an inhibitor of the protein-protein interaction between murine double minute 2 (MDM2) and p53.<sup>1</sup> It induces apoptosis in WaGa and PeTa Merkel cell carcinoma cells when used at a concentration of 100 nM. DS-3032 (1 μM) decreases the viability of p53 wild-type MKL-1 cells but not p53 knockout MLK-1 cells. *In vivo*, DS-3032 (100 mg/kg per day) reduces tumor volume and increases survival in a patient-derived xenograft (PDX) mouse model of Merkel cell carcinoma.

## Reference

1. Ananthapadmanabhan, V., Frost, T.C., Soroko, K.M., *et al.* Milademetan is a highly potent MDM2 inhibitor in Merkel cell carcinoma. *JCI Insight* **7**(13), e160513 (2022).

### 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, 10/09/2024

## 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)