

PRODUCT INFORMATION



SJ 172550

Item No. 19999

CAS Registry No.: 431979-47-4
Formal Name: 2-[2-chloro-4-[(1,5-dihydro-3-methyl-5-oxo-1-phenyl-4H-pyrazol-4-ylidene)methyl]-6-ethoxyphenoxy]-acetic acid, methyl ester

Synonym: MDMX Inhibitor II

MF: C₂₂H₂₁ClN₂O₅

FW: 428.9

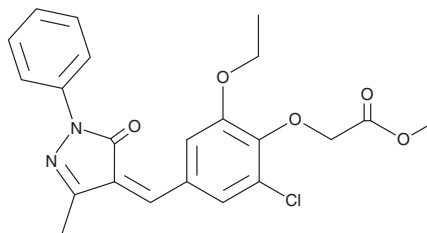
Purity: ≥98%

UV/Vis.: λ_{max}: 251, 343 nm

Supplied as: A crystalline 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

SJ 172550 is supplied as a crystalline solid. A stock solution may be made by dissolving the SJ 172550 in the solvent of choice. SJ 172550 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas, and slightly soluble in ethanol. The solubility of SJ 172550 in DMSO and DMF is approximately 10 and 30 mg/ml, respectively.

SJ 172550 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SJ 172550 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SJ 172550 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SJ 172550 is an inhibitor of MDMX that disrupts MDMX-p53 peptide interaction with an EC₅₀ value of 4.3 μM.¹ It forms a covalent, but reversible, cysteine adduct in the MDMX p53-binding domain, locking MDMX into a conformation that prevents binding to p53.² SJ 172550 induces apoptosis, increases p53 target gene expression, and causes p53-dependent cell death in retinoblastoma cells.¹

References

1. Reed, D.J., Shen, Y., Shelat, A.A., *et al.* Identification and characterization of the first small molecule inhibitor of MDMX. *J. Biol. Chem.* **285**(4), 10786-10796 (2010).
2. Bista, M., Smithson, D., Pecak, A., *et al.* On the mechanism of action of SJ-172550 in inhibiting the interaction of MDM4 and p53. *PLoS One* **7**(6), e37518 (2012).

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