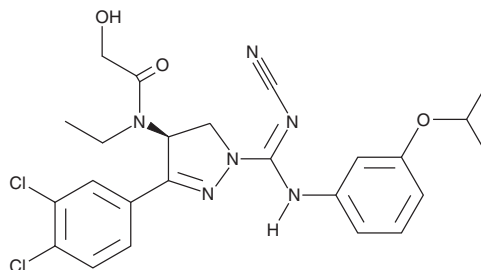


PRODUCT INFORMATION



(S)-BAY-598 Item No. 18238

CAS Registry No.: 1906919-67-2
Formal Name: (S,Z)-N-(1-(N'-cyano-N-(3-(difluoromethoxy)phenyl)carbamidoyl)-3-(3,4-dichlorophenyl)-4,5-dihydro-1H-pyrazol-4-yl)-N-ethyl-2-hydroxyacetamide
MF: C₂₂H₂₀Cl₂F₂N₆O₃
FW: 525.3
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 324 nm
Supplied as: A crystalline solid
Storage: 2-8°C
Stability: ≥4 years



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

Laboratory Procedures

(S)-BAY-598 is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-BAY-598 in the solvent of choice. (S)-BAY-598 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of (S)-BAY-598 in these solvents is approximately 30 mg/ml.

Description

(S)-BAY-598 is a selective inhibitor of SMYD2. It inhibits the methylation of p53 at lysine 370 *in vitro* and in whole cells with IC₅₀ values of 27 nM and <1 μM, respectively.¹ (S)-BAY-598 exhibits 100-fold selectivity for SMYD2 over other histone methyltransferases and non-epigenetic enzymes. For more information on (S)-BAY-598 please visit the Structural Genomics Consortium (SGC). The negative control, BAY-369, for (S)-BAY-598 is also available exclusively through the SGC.

Reference

1. Eggert, E., Hillig, R.C., Koehr, S., *et al.* Discovery and characterization of a highly potent and selective aminopyrazoline-based *in vivo* probe (BAY-598) for the protein lysine methyltransferase SMYD2. *J. Med. Chem.* **59**(10), 4578-4600 (2016).

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, 11/21/2022

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