

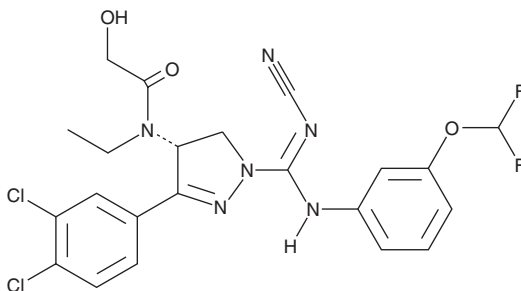
PRODUCT INFORMATION



(R)-BAY-598

Item No. 29174

CAS Registry No.: 1906920-28-2
Formal Name: N-[(4R)-1-[(cyanoamino) [[3-(difluoromethoxy)phenyl]imino] methyl]-3-(3,4-dichlorophenyl)-4,5-dihydro-1H-pyrazol-4-yl]-N-ethyl-2-hydroxy-acetamide
MF: C₂₂H₂₀Cl₂F₂N₆O₃
FW: 525.3
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

(R)-BAY-598 is supplied as a solid. A stock solution may be made by dissolving the (R)-BAY-598 in the solvent of choice, which should be purged with an inert gas. (R)-BAY-598 is soluble in the organic solvent DMSO.

Description

(R)-BAY-598 is an inhibitor of the lysine N-methyltransferase SMYD2 (IC₅₀ = 1.7 μM) and a less active enantiomer of (S)-BAY-598 (Item No. 18238).¹ It is selective for SMYD2 over proteinase-activated receptor 1 (PAR1; IC₅₀ = >30 μM).

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, 12/01/2022

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