

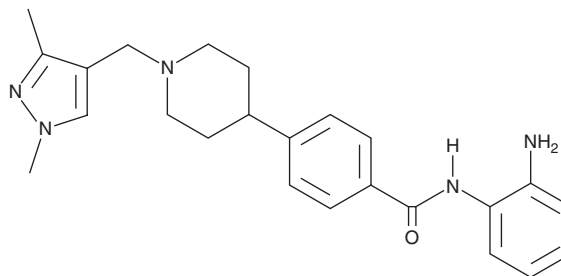
PRODUCT INFORMATION



CXD101

Item No. 21159

CAS Registry No.: 934828-12-3
Formal Name: N-(2-aminophenyl)-4-[1-[(1,3-dimethyl-1H-pyrazol-4-yl)methyl]-4-piperidinyl]-benzamide
Synonym: AZD 9468
MF: C₂₄H₂₉N₅O
FW: 403.5
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 298 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

CXD101 is supplied as a crystalline solid. A stock solution may be made by dissolving the CXD101 in the solvent of choice, which should be purged with an inert gas. CXD101 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CXD101 in ethanol is approximately 1 mg/ml and approximately 10 mg/ml in DMSO and DMF.

CXD101 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CXD101 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CXD101 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CXD101 is a histone deacetylase (HDAC) inhibitor with structural similarities to mocetinostat (Item No. 18287), MS-275 (Item No. 13284), and JNJ-26481585 (Item No. 14088). The biological activity of CXD101 has not been published.

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