

# PRODUCT INFORMATION



## CCS-1477

Item No. 34546

**CAS Registry No.:** 2222941-37-7  
**Formal Name:** (6S)-1-(3,4-difluorophenyl)-6-[5-(3,5-dimethyl-4-isoxazolyl)-1-(trans-4-methoxycyclohexyl)-1H-benzimidazol-2-yl]-2-piperidinone

**Synonym:** CBP-IN-1

**MF:** C<sub>30</sub>H<sub>32</sub>F<sub>2</sub>N<sub>4</sub>O<sub>3</sub>

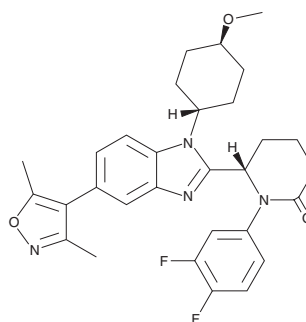
**FW:** 534.6

**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

CCS-1477 is supplied as a solid. A stock solution may be made by dissolving the CCS-1477 in the solvent of choice, which should be purged with an inert gas. CCS-1477 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CCS-1477 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

CCS-1477 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CCS-1477 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CCS-1477 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

CCS-1477 is an inhibitor of p300 and CREB-binding protein (CBP) histone acetyltransferases.<sup>1</sup> It selectively binds to p300 and CBP over bromodomain-containing protein 4 (BRD4; K<sub>d</sub>s = 1.3, 1.7, and 222 nM, respectively). CCS-1477 inhibits the proliferation of OPM-2 multiple myeloma cells (GI<sub>50</sub> = 5 nM). It induces tumor regression in MOLM-16 acute myeloid leukemia (AML) and OPM-2 mouse xenograft models when administered at a dose of 20 mg/kg.

## Reference

1. Brooks, N., Raja, M., Young, B.W., *et al.* CCS1477: A novel small molecule inhibitor of p300/CBP bromodomain for the treatment of acute myeloid leukaemia and multiple myeloma. *Blood* 134(Supplement 1), 2560 (2019).

### 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, 12/12/2022

## 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