

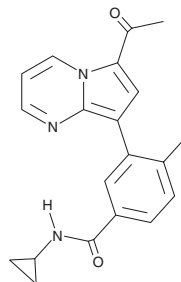
# PRODUCT INFORMATION



TP-472

Item No. 20030

**CAS Registry No.:** 2079895-62-6  
**Formal Name:** 3-(6-acetylpyrrolo[1,2-a]pyrimidin-8-yl)-N-cyclopropyl-4-methylbenzamide  
**MF:** C<sub>20</sub>H<sub>19</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 333.4  
**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

TP-472 is supplied as a solid. A stock solution may be made by dissolving the TP-472 in the solvent of choice, which should be purged with an inert gas. TP-472 is soluble in the organic solvent DMSO.

## Description

TP-472 is an inhibitor of bromodomain BRD9 ( $K_d = 33$  nM;  $EC_{50} = 320$  nM in a NanoBRET assay). It has >30-fold selectivity over all other bromodomain family members with the exception of the highly homologous BRD7 bromodomain ( $K_d = 340$  nM in an isothermal titration calorimetry (ITC) assay). See the Structural Genomic Consortium (SGC) website for more information.

### 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, 11/30/2022

## 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