

# PRODUCT INFORMATION



**AU-15330**

Item No. 43318

**CAS Registry No.:** 2380274-50-8

**Formal Name:** (2S,4R)-1-((S)-2-(2-(4-(3-amino-6-(2-hydroxyphenyl)pyridazin-4-yl)piperazin-1-yl)acetamido)-3,3-dimethylbutanoyl)-4-hydroxy-N-((S)-1-(4-(4-methylthiazol-5-yl)phenyl)ethyl)pyrrolidine-2-carboxamide

**MF:** C<sub>39</sub>H<sub>49</sub>N<sub>9</sub>O<sub>5</sub>S

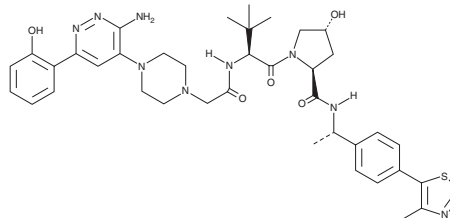
**FW:** 755.9

**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

AU-15330 is supplied as a solid. A stock solution may be made by dissolving the AU-15330 in the solvent of choice, which should be purged with an inert gas. AU-15330 is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

## Description

AU-15330 is a proteolysis-targeting chimera (PROTAC) that contains a von Hippel-Lindau (VHL) ligand conjugated to AU-15139, a ligand of the Brahma homolog (BHM) and Brahma-related gene 1 (BRG1), also known as SMARCA2 and SMARCA4, respectively.<sup>1</sup> It induces degradation of BRM and BRG1 in HeLa cervical cancer cells when used at a concentration of 1 μM but also induces degradation of polybromo-1D (PBRM1) at the same concentration. AU-15330 decreases proliferation in a panel of 14 cancer cell lines (IC<sub>50</sub>s = 0.392-28.85 nM). It reduces protein levels of forkhead box protein A1 (FOXA1), ETS-related gene (ERG), Myc, and prostate-specific antigen in VCaP prostate cancer cells when used at a concentration of 1 μM for 24 hours. AU-15330 reduces tumor volume in a VCaP mouse xenograft model.

## Reference

1. Xiao, L., Parolia, A., Qiao, Y., *et al.* Targeting SWI/SNF ATPases in enhancer-addicted prostate cancer. *Nature* **601**(7893), 434-439 (2022).

### 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, 05/02/2025

## 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](http://WWW.CAYMANCHEM.COM)