

# PRODUCT INFORMATION

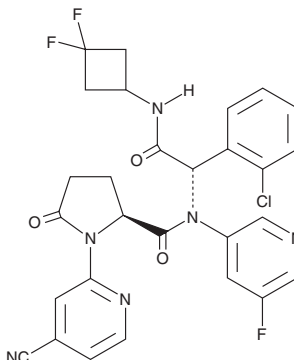


## AG-120

Item No. 19894

**CAS Registry No.:** 1448347-49-6  
**Formal Name:** (2S)-1-(4-cyano-2-pyridinyl)-5-oxo-L-prolyl-2-(2-chlorophenyl)-N-(3,3-difluorocyclobutyl)-N<sup>2</sup>-(5-fluoro-3-pyridinyl)-glycinamide

**Synonym:** Ivosidenib  
**MF:** C<sub>28</sub>H<sub>22</sub>ClF<sub>3</sub>N<sub>6</sub>O<sub>3</sub>  
**FW:** 583.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 246, 307 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

AG-120 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-120 in the solvent of choice. AG-120 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of AG-120 in these solvents is approximately 30 mg/ml.

### Description

AG-120 is an orally available inhibitor of isocitrate dehydrogenase type 1 (IDH1) that specifically inhibits a mutated form of IDH1 in the cytoplasm, preventing the formation of 2-hydroxyglutarate.<sup>1</sup> AG-120 may induce cellular differentiation as well as inhibit proliferation in IDH1-expressing tumor cells.<sup>1</sup>

### Reference

1. Gu, C.-H. Pharmaceutical compositions of therapeutically active compounds. World Intellectual Property Organization - International Bureau Publication WO 2015/138839 A1 (2015), 1-13, PCT/US2015/020349.

#### 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, 09/01/2023

#### 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)