

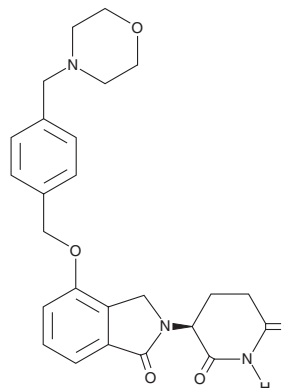
PRODUCT INFORMATION



Iberdomide

Item No. 22615

CAS Registry No.: 1323403-33-3
Formal Name: 3S-[1,3-dihydro-4-[[4-(4-morpholinylmethyl)phenyl]methoxy]-1-oxo-2H-isoindol-2-yl]-2,6-piperidinedione
Synonym: CC-220
MF: C₂₅H₂₇N₃O₅
FW: 449.5
Purity: ≥98%
UV/Vis.: λ_{max}: 284 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

Iberdomide is supplied as a crystalline solid. A stock solution may be made by dissolving the iberdomide in the solvent of choice, which should be purged with an inert gas. Iberdomide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of iberdomide in these solvents is approximately 12 mg/ml.

Iberdomide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, iberdomide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Iberdomide has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Iberdomide is a modulator of cereblon, a substrate receptor in the CRL4 ubiquitin ligase complex, that has an IC₅₀ value of 60 nM in a competitive TR-FRET assay.¹ In the same assay, it has a higher affinity for cereblon than the related modulators lenalidomide (Item No. 14643; IC₅₀ = 1,500 nM) and pomalidomide (Item No. 19877; IC₅₀ = 1,200 nM). Iberdomide also has a greater potency for degrading the CRL4 complex substrates Ikaros and Aiolos (EC₅₀s = 1 and 0.5 nM) but has no degradative activity for GSPT1 or CK1α.

Reference

1. Matyskiela, M.E., Zhang, W., Man, H.W., *et al.* A cereblon modulator (CC-220) with improved degradation of Ikaros and Aiolos. *J. Med. Chem.* (2017).

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, 11/17/2022

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