

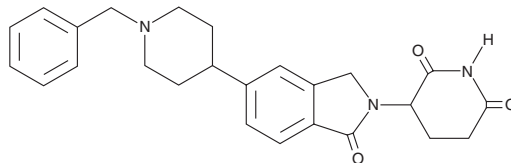
# PRODUCT INFORMATION



## NVP-DKY709

Item No. 41295

CAS Registry No.: 2291360-73-9  
Formal Name: 3-[1,3-dihydro-1-oxo-5-[1-(phenylmethyl)-4-piperidinyl]-2H-isoindol-2-yl]-2,6-piperidinedione  
MF:  $C_{25}H_{27}N_3O_3$   
FW: 417.5  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

NVP-DKY709 is supplied as a solid. A stock solution may be made by dissolving the NVP-DKY709 in the solvent of choice, which should be purged with an inert gas. NVP-DKY709 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

NVP-DKY709 is a degrader of Ikaros family zinc finger protein 2 (IKZF2) with a half-maximal degradation concentration ( $DC_{50}$ ) value of 4 nM.<sup>1</sup> It is selective for IKZF2 over 8,656 other proteins, including IKZF1, IKZF3, and IKZF5, at 2.5  $\mu\text{M}$ , but it does degrade IKZF4, which has the same zinc finger 2  $\beta$ -hairpin sequence as IKZF2, and Sal-like protein 4 (SALL4;  $DC_{50}$ s = 13 and 2 nM, respectively). It increases IL-2 secretion induced by phytohemagglutinin in Jurkat T cells in a concentration-dependent manner. NVP-DKY709 (100 mg/kg), alone or in combination with an antibody against programmed death protein 1 (PD-1), decreases tumor volume in an MDA-MB-231 humanized mouse xenograft model.

### Reference

1. Bonazzi, S., d'Hennezel, E., Beckwith, R.E., *et al.* Discovery and characterization of a selective IKZF2 glue degrader for cancer immunotherapy. *Cell Chem. Biol.* **30(3)**, 235-247.e12 (2023).

#### 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, 09/5/2024

#### 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