

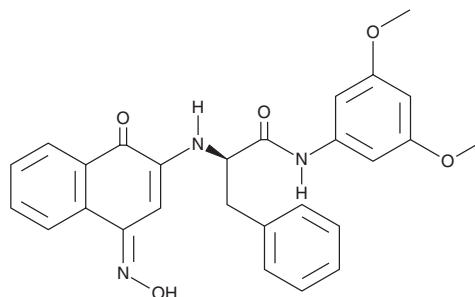
# PRODUCT INFORMATION



**CAY10763**

Item No. 30593

**CAS Registry No.:** 2364458-49-9  
**Formal Name:** (αR)-α-[[1,4-dihydro-4-(hydroxyimino)-1-oxo-2-naphthalenyl]amino]-N-(3,5-dimethoxyphenyl)-benzenepropanamide  
**MF:** C<sub>27</sub>H<sub>25</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 471.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 242 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

## Laboratory Procedures

CAY10763 is supplied as a solid. A stock solution may be made by dissolving the CAY10763 in the solvent of choice, which should be purged with an inert gas. CAY10763 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CAY10763 in these solvents is approximately 30 mg/ml.

## Description

CAY10763 is a dual inhibitor of indoleamine 2,3-dioxygenase 1 (IDO1; IC<sub>50</sub> = 46 nM) and STAT3 activation.<sup>1</sup> It binds to STAT3 (K<sub>d</sub> = 530 nM) and selectively reduces the levels of STAT3 phosphorylated at the tyrosine in position 705 (STAT3<sup>Y705</sup>) over phosphorylated STAT3<sup>S727</sup>, STAT1, and STAT5 in SKOV3 cells when used at a concentration of 500 nM. It also inhibits STAT3 nuclear translocation in SKOV3 cells. CAY10763 is cytotoxic to HCT116, SKOV3, A549, and HepG2 cancer cells (IC<sub>50</sub>s = 37, 28, 33, and 12 nM, respectively). It reduces tumor growth in a B16/F10 murine melanoma model and a HepG2 mouse xenograft model when administered at doses of 100 and 10 mg/kg, respectively.

## Reference

1. Huang, R., Jing, X., Huang, X., *et al.* Bifunctional naphthoquinone aromatic amide-oxime derivatives exert combined immunotherapeutic and antitumor effects through simultaneous targeting of indoleamine-2,3-dioxygenase and signal transducer and activator of transcription 3. *J. Med. Chem.* **63**(4), 1544-1563 (2020).

### 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, 01/15/2021

## 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