

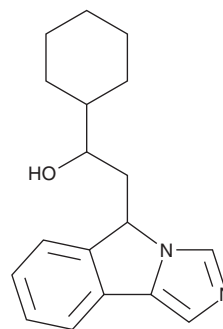
PRODUCT INFORMATION



NLG-919 analog

Item No. 21509

CAS Registry No.: 1402836-58-1
Formal Name: α -cyclohexyl-5H-imidazo[5,1-a]isoindole-5-ethanol
MF: C₁₈H₂₂N₂O
FW: 282.4
Purity: \geq 98% (mixture of enantiomers)
UV/Vis.: λ_{max} : 213, 274 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

NLG-919 analog is supplied as a crystalline solid. A stock solution may be made by dissolving the NLG-919 analog in the solvent of choice, which should be purged with an inert gas. NLG-919 analog is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NLG-919 analog in ethanol is approximately 14 mg/ml and approximately 16 mg/ml in DMSO and DMF.

NLG-919 analog is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NLG-919 analog should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NLG-919 analog has a solubility of approximately 0.3 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NLG-919 analog is an inhibitor of indoleamine 2,3-dioxygenase 1 (IDO1) and tryptophan 2,3-dioxygenase (TDO; IC₅₀s = 79 and 25 nM, respectively, in cell-free assays).¹ It decreases kynurenine levels in cells expressing human IDO1 (EC₅₀ = 83 nM).² NLG-919 analog (0.025-1 μ M) decreases the percentage of anti-CD3- and anti-CD28-activated Jurkat T cells that are positive for programmed cell death protein 1 (PD-1).³ *In vivo*, NLG-919 analog (100 mg/kg) decreases tumor volume and weight and the percentage of intratumoral CD4⁺CD25⁺ regulatory T cells (Tregs) in a B16/F10 murine melanoma model when administered alone or in combination with paclitaxel (Item No. 10461).⁴

Reference

1. Seegers, N., van Doornmalen, A.M., Uitdehaag, J.C., *et al.* High-throughput fluorescence-based screening assays for tryptophan-catabolizing enzymes. *J. Biomol. Screen.* **19(9)**, 1266-1274 (2014).
2. Kumar, S., Waldo, J.P., Jaipuri, F.A., *et al.* Discovery of clinical candidate (1R,4r)-4-((R)-2-((S)-6-fluoro-5H-imidazo[5,1-a]isoindol-5-yl)-1-hydroxyethyl)cyclohexan-1-ol (navoximod), a potent and selective inhibitor of indoleamine 2,3-dioxygenase 1. *J. Med. Chem.* **62(14)**, 6705-3766 (2019).
3. Qin, R., Zhao, C., Wang, C.J., *et al.* Tryptophan potentiates CD8⁺ T cells against cancer cells by TRIP12 tryptophanylation and surface PD-1 downregulation. *J. Immunother. Cancer* **9(7)**, e002840 (2021).
4. Meng, X., Du, G., Ye, L., *et al.* Combinatorial antitumor effects of indoleamine 2,3-dioxygenase inhibitor NLG919 and paclitaxel in a murine B16-F10 melanoma model. *Int. J. Immunopathol. Pharmacol.* **30(3)**, 215-226 (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, 10/23/2024

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