

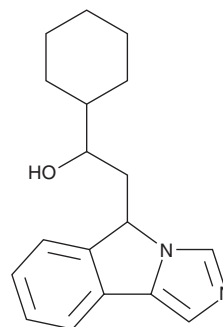
PRODUCT INFORMATION



(±)-NLG-919

Item No. 21509

CAS Registry No.: 1402836-58-1
Formal Name: α-cyclohexyl-5H-imidazo[5,1-a]isoindole-5-ethanol
Synonym: GDC-0919
MF: C₁₈H₂₂N₂O
FW: 282.4
Purity: ≥98% (mixture of enantiomers)
UV/Vis.: λ_{max}: 213, 274 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

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

(±)-NLG-919 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (±)-NLG-919 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (±)-NLG-919 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 is an indoleamine 2,3-dioxygenase (IDO) pathway inhibitor ($K_i = 7$ nM; $EC_{50} = 75$ nM).¹ (±)-NLG-919 inhibits IDO-induced T cell suppression and restores T cell response during allogeneic mixed lymphocyte reactions *in vitro* (ED_{50} s = 80 and 120 nM, respectively). *In vivo*, (±)-NLG-919 induces a 95% reduction in tumor volume when administered in conjunction with a pmel-1 T cell vaccine.

Reference

1. Mautino, M.R., Jaipuri, F.A., Waldo, J., *et al.* Abstract 491: NLG919, a novel indoleamine-2,3-dioxygenase (IDO)-pathway inhibitor drug candidate for cancer therapy. *Cancer Res.* **73**(8 Suppl.), 491 (2013).

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, 11/07/2022

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