# **PRODUCT** INFORMATION V-125



′532

CAS Registry No.:	1807740-94-8		
Formal Name:	6-[ethyl(5,6,7,8-tetrahydro-3,5,5,8,8-		
	pentamethyl-2-naphthalenyl)aminol-		
	3-pyridinecarboxylic acid		
MF:	$C_{23}H_{30}N_2O_2$		
FW:	366.5		
Purity:	≥98%		
UV/Vis.:	λ <sub>max</sub> : 291 nm	$X \sim \cdot \sim H$	
Supplied as:	A solid	/ \ 0	
Storage:	-20°C		
Stability:	≥2 years		
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.			

### Laboratory Procedures

V-125 is supplied as a solid. A stock solution may be made by dissolving the V-125 in the solvent of choice, which should be purged with an inert gas. V-125 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). V-125 is soluble in DMF at a concentration of approximately 2 mg/ml. V-125 is slightly soluble in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of V-125 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of V-125 in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

V-125 is an agonist of retinoid X receptors (RXRs).<sup>1</sup> It induces homodimerization of RXR in HCT116 cells in a two-hybrid assay (EC<sub>50</sub> = 13.8 nM). V-125 selectively induces transactivation of RXRa in human RXRα-overexpressing HCT116 cells over transactivation of RARα in human RARα-overexpressing HEK293 cells when used at concentrations of 25 and 100 nM, respectively. In vivo, V-125 (80 mg/kg in the diet) inhibits vinyl carbamate-induced lung tumor growth in mice.<sup>2</sup> Unlike the RXR agonist bexarotene (Item No. 11571), V-125 does not increase plasma triglyceride or cholesterol levels in mice.

### References

- 1. Heck, M.C., C.E., W., Shahani, P.H., et al. Modeling, synthesis, and biological evaluation of potential retinoid X receptor (RXR)-selective agonists: Analogues of 4 [1-(3,5,5,8,8-pentamethyl-5,6,7,8-tetrahydro-2-naphthyl)ethynyl]benzoic acid (bexarotene) and 6 (ethyl(5,5,8,8-tetrahydronaphthalen-2-yl)amino) nicotinic acid (NEt-TMN). J. Med. Chem. 59(19), 8924-8940 (2016).
- 2. Reich, L.A., Moerland, J.A., Leal, A.S., et al. The rexinoid V 125 reduces tumor growth in preclinical models of breast and lung cancer. Sci. Rep. 12(1), 293 (2022).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

### SAFFTY 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

uyer 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, 11/11/2022

## 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