# **PRODUCT** INFORMATION



Narciclasine

Item No. 20361

CAS Registry No.:	29477-83-6	
Formal Name:	(2S,3R,4S,4aR)-3,4,4a,5-tetrahydro-2,3,4,7-	ŎН
	tetrahydroxy-[1,3]dioxolo[4,5-j]phenanthridin-6(2H)-one	
Synonyms:	(+)-Lycoricidinol, (+)-Narciclasine, NSC 266535	
MF:	C <sub>14</sub> H <sub>13</sub> NO <sub>7</sub>	н
FW:	307.3	
Purity:	≥98%	
UV/Vis.:	λ <sub>max</sub> : 213, 252, 302 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	
Stability:	≥4 years	OH
Item Origin:	Plant/Narcissus sp.	

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

# Laboratory Procedures

Narciclasine is supplied as a crystalline solid. A stock solution may be made by dissolving the narciclasine in the solvent of choice, which should be purged with an inert gas. Narciclasine is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

# Description

Narciclasine is an isocarbostyril alkaloid that has been found in Narcissus and has diverse biological activities.<sup>1-5</sup> It selectively induces apoptosis in PC3 prostate and MCF-7 and MDA-MB-231 breast cancer cells over CCD-25Lu fibroblasts at 1  $\mu$ M.<sup>1</sup> Narciclasine has antiproliferative activity against a panel of six glioblastoma cancer cell lines (mean  $IC_{50}$  = ~50 nM), induces actin stress fiber formation in U373 MG glioblastoma cells, and increases survival in an orthotopic Hs 683 glioblastoma mouse xenograft model.<sup>2</sup> It also inhibits neurite outgrowth in SH-SY5Y cells with a 10% effective concentration (EC<sub>10</sub>) value of 4.44 nM and suppresses the proliferation of isolated human T cells ( $IC_{50} = 14 \text{ nM}$ ).<sup>3,4</sup> Intradermal administration of narciclasine (2 mg/kg) decreases disease severity and prevents increases in spleen weight in a mouse model of psoriasis induced by imiquimod (Item No. 14956).<sup>5</sup>

# Reference

- 1. Dumont, P., Ingrassia, L., Rouzeau, S., et al. The Amaryllidaceae isocarbostyril narciclasine induces apoptosis by activation of the death receptor and/or mitochondrial pathways in cancer cells but not in normal fibroblasts. Neoplasia 9(9), 766-776 (2007).
- 2. Lefranc, F., Sauvage, S., Van Goietsenoven, G., et al. Narciclasine, a plant growth modulator, activates Rho and stress fibers in glioblastoma cells. Mol. Cancer Ther. 8(7), 1739-1750 (2009).
- 3. Braun, G., Herberth, G., Krauss, M., et al. Neurotoxic mixture effects of chemicals extracted from blood of pregnant women. Science 386(6719), 301-309 (2024).
- 4. Wang, W.-L., Wu, X.-Y., Luo, X.-Y., et al. Immunosuppressive alkaloids from Narcissus tazetta subsp. Chinensis and the mechanism of (+)-narciclasine in vitro and in vivo. Phytochemistry 225:114198, (2024).
- 5. Kong, Y., Jiang, J., Huang, Y., et al. Narciclasine inhibits phospholipase A2 and regulates phospholipid metabolism to ameliorate psoriasis-like dermatitis. Front. Immunol. 13:1094375, (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

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, 12/12/2024

# 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