# **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-             | UH :    |
|                   | tetrahydroxy-[1,3]dioxolo[4,5-j]phenanthridin-6(2H)-one | OH      |
| Synonyms:         | (+)-Lycoricidinol, (+)-Narciclasine                     | ÍI Í    |
| MF:               | C <sub>14</sub> H <sub>13</sub> NO <sub>7</sub>         |         |
| FW:               | 307.3   |         |
| Purity:           | ≥98%  |         |
| UV/Vis.:          | λ <sub>max</sub> : 213, 252, 302 nm                     | N N     |
| Supplied as:      | A crystalline solid                                     | т Т П н |
| Storage:          | -20°C   | I O     |
| Stability:        | ≥4 years  |         |
|                   |   |         |

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. Narciclasine is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 100 mM.

### Description

Narciclasine is a plant growth inhibitor that can be isolated from Narcissus bulbs. At 1 µM, it has been shown to induce apoptosis-mediated cytotoxicity in human cancer cells in vitro but not in normal fibroblasts.<sup>1</sup> Narciclasine has been shown to regulate the Rho/Rho kinase/LIM kinase/cofilin signaling pathway by increasing GTPase RhoA activity, as well as inducing actin stress fiber formation in a RhoA-dependent manner.<sup>1</sup>

#### Reference

1. 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).

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.

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, 10/31/2022

## 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