# **PRODUCT** INFORMATION



## Gedunin

Item No. 17328

CAS Registry No.:	2753-30-2	0
Formal Name:	(1S,3aS,4aR,4bS,5R,6aR,10aR,10bR,12aS)-5-	
	(acetyloxy)-1-(3-furanyl)-1,5,6,6a,7,10a,10b,11,12,12a-	
	decahydro-4b,7,7,10a,12a-pentamethyl-oxireno[c]	$\checkmark$
	phenanthro[1,2-d]pyran-3,8(3aH,4bH)-dione	
Synonym:	NSC 113497	o o
MF:	$C_{28}H_{34}O_7$	
FW:	482.6	
Purity:	≥98%	Н О
UV/Vis.:	λ <sub>max</sub> : 220 nm	
Supplied as:	A crystalline solid	$\downarrow$ $\downarrow$ $\downarrow$ $\downarrow$ $\downarrow$ $\downarrow$
Storage:	-20°C	
Stability:	≥2 years	0^

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

#### Laboratory Procedures

Gedunin is supplied as a crystalline solid. A stock solution may be made by dissolving the gedunin in the solvent of choice. Gedunin is soluble in organic solvents such as ethanol, methanol, and acetonitrile, which should be purged with an inert gas. The solubility of gedunin in these solvents is approximately 1 mg/ml.

#### Description

Gedunin is a natural inhibitor of the heat shock protein (HSP90; Item Nos. 22734 | 22735) co-chaperone p23 that also inhibits HSP90 expression in human teratocarcinomal NTERA-2 cells in vitro at 5  $\mu$ g/ml.<sup>1</sup> This tetranortriterpenoid, which is isolated from A. indica, binds to and blocks the chaperone activity of p23 to induce apoptosis in HeLa-PR<sub>B</sub> cells in vitro at a concentration of 20  $\mu$ M.<sup>2</sup> Gedunin inhibits breast cancer cell proliferation in vitro, with IC50 values of 8.84 and 3.22 µM in MCF-7 and SKBr-3 cells, respectively, and inhibits the growth of PANC-1 pancreatic cancer cells (IC<sub>50</sub> = 25  $\mu$ M) by targeting the sonic hedgehog pathway to induce apoptosis.<sup>3,4</sup> It also exerts anti-inflammatory effects in vivo. In a mouse model of articular inflammation induced by zymosan (Item No. 21175), gedunin (0.05-0.5 mg/kg, i.p.) reduces edema formation and the production of inflammatory cytokines.<sup>5</sup> Gedunin (0.5 mg/kg) also inhibits the pleural accumulation of eosinophils and activated T lymphocytes in an ovalbumin-sensitized mouse model of allergic inflammation when administered prior to ovalbumin rechallenge.<sup>6</sup> It also targets the lipopolysaccharide binding site and thus blocks Toll-like receptor 4 (TLR4) signaling in macrophages in vitro at a concentration of 10  $\mu$ M.<sup>7</sup>

#### References

- 1. Tharmarajah, L., Samarakoon, S.R., Ediriweera, M.K., et al. Biomed. Res. Int. 2017, 2413197 (2017).
- 2. Patwardhan, C.A., Fauq, A., Peterson, L.B., et al. J. Biol. Chem. 288(10), 7313-7325 (2013).
- 3. Brandt, G.E.L., Schmidt, M.D., Prisinzano, T.E., et al. J. Med. Chem. 51(20), 6495-6502 (2008).
- 4. Subramani, R., Gonzalez, E., Nandy, S.B., et al. Oncotarget 8(7), 10891-10904 (2017).
- 5. Conte, F.P., Kerraris, F.K., Costa, T.E.M.M., et al. Molecules 20(2), 2636-2657 (2015).
- 6. Ferraris, F.K., Moret, K.H., Figueiredo, A.B.C., et al. Int. Immunopharmacol. 14(1), 82-93 (2012).
- 7. Borges, P.V., Moret, K.H., Maya-Monteiro, C.M., et al. Mol. Pharm. 88(5), 949-961

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

#### SAFFTY DATA

al 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, 01/18/2018

### 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