

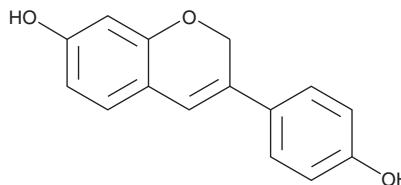
# PRODUCT INFORMATION



## Idronoxil

Item No. 30783

CAS Registry No.: 81267-65-4  
Formal Name: 3-(4-hydroxyphenyl)-2H-1-benzopyran-7-ol  
Synonyms: Haginin E, Phenoxodiol  
MF:  $C_{15}H_{12}O_3$   
FW: 240.3  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: Synthetic



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

### Laboratory Procedures

Idronoxil is supplied as a solid. A stock solution may be made by dissolving the idronoxil in the solvent of choice, which should be purged with an inert gas. Idronoxil is soluble in the organic solvent DMSO.

### Description

Idronoxil is a phenol and derivative of genistein (Item No. 10005167) that has been found in *B. tournefortii* and has anticancer activity.<sup>1-3</sup> It reduces the viability of R182S, R127, Hey, CP70, A2780, R187, R188, and R207 primary ovarian cancer cells, but not non-cancerous ovarian surface epithelial (OSE) cells, when used at a concentration of 10  $\mu\text{g}/\text{ml}$ .<sup>2</sup> Idronoxil (1  $\mu\text{g}/\text{ml}$ ) reduces colony formation and induces apoptosis in R127 and CP70 primary ovarian cancer cells, respectively, and restores sensitivity to Fas-mediated apoptosis in CP70 cells. *In vivo*, idronoxil (50 and 75 mg/kg) increases latency to tumor formation and reduces tumor multiplicity in a rat model of mammary carcinogenesis induced by dimethylbenz[a]anthracene (DMBA).<sup>3</sup>

### References

1. Rahmani, R., Bouajila, J., Jouaidi, M., *et al.* African mustard (*Brassica tournefortii*) as source of nutrients and nutraceuticals properties. *J. Food Sci.* **85**(6), 1856-1871 (2020).
2. Kamsteeg, M., Rutherford, T., Sapi, E., *et al.* Phenoxodiol – an isoflavone analog – induces apoptosis in chemoresistant ovarian cancer cells. *Oncogene* **22**(17), 2611-2620 (2003).
3. Constantinou, A.I., Mehta, R., and Husband, A. Phenoxodiol, a novel isoflavone derivative, inhibits dimethylbenz[a]anthracene (DMBA)-induced mammary carcinogenesis in female Sprague-Dawley rats. *Eur. J. Cancer* **39**(7), 1012-1018 (2004).

#### 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, 12/13/2022

#### 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