

PRODUCT INFORMATION



Acetylharpagide

Item No. 39930

CAS Registry No.: 6926-14-3

Formal Name: (1S,4aS,5R,7S,7aS)-7-(acetyloxy)-1,4a,5,6,7,7a-hexahydro-4a,5-dihydroxy-7-methylcyclopenta[c]pyran-1-yl, β -D-glucopyranoside

Synonyms: 8-Acetylharpagide, 8-O-Acetylharpagide, 8-AcHarp, Harpagide Acetate

MF: $C_{17}H_{26}O_{11}$

FW: 406.4

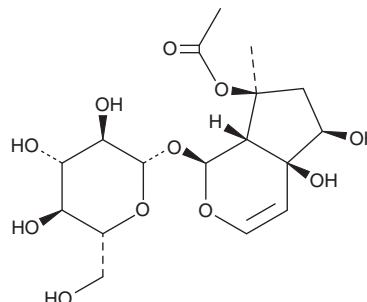
Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Scrophularia ningpoensis* Hemsl.



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

Laboratory Procedures

Acetylharpagide is supplied as a solid. A stock solution may be made by dissolving the acetylharpagide in the solvent of choice, which should be purged with an inert gas. Acetylharpagide is soluble in methanol. Acetylharpagide is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Acetylharpagide is an iridoid glycoside that has been found in *A. decumbens* and has diverse biological activities.¹⁻³ It reduces the proliferation of A431 skin cancer cells ($IC_{50} = 310 \mu\text{M}$).¹ Acetylharpagide (0.005% in the drinking water) reduces the number of tumors per mouse and the percentage of mice with tumors in a mouse model of 4-nitroquinolin 1-oxide-induced carcinogenesis.² It decreases carrageenan-induced paw swelling in rats when administered at a dose of 10 mg/kg.³ Acetylharpagide (10 mg/kg) reduces blood glucose levels in a mouse model of type I diabetes induced by alloxan (Item No. 9002196).

References

1. Kuria, K.A.M., Chepkwony, H., Govaerts, C., *et al.* The antiplasmodial activity of isolates from *Ajuga remota*. *J. Nat. Prod.* **65**(5), 789-793 (2002).
2. Takasaki, M., Tokuda, H., Nishino, H., *et al.* Cancer chemopreventive agents (antitumor-promoters) from *Ajuga decumbens*. *J. Nat. Prod.* **62**(7), 972-975 (1999).
3. Ahmed, B., Al-Rehaily, A.J., Al-Howiriny, T.A., *et al.* Scropolioside-D₂ and harpagoside-B: Two new iridoid glycosides from *Scrophularia deserti* and their antidiabetic and antiinflammatory activity. *Biol. Pharm. Bull.* **26**(4), 462-467 (2003).

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, 02/29/2024

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