

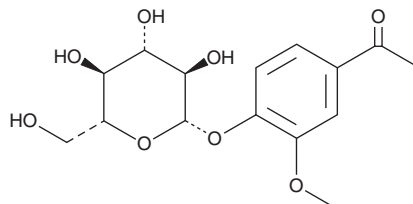
PRODUCT INFORMATION



Androsin

Item No. 36191

CAS Registry No.: 531-28-2
Formal Name: 1-[4-(β-D-glucopyranosyloxy)-3-methoxyphenyl]-ethanone
MF: C₁₅H₂₀O₈
FW: 328.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 268, 302 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Nerium indicum*



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

Laboratory Procedures

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

Description

Androsin is a phenol glucoside that has been found in *P. kurroa* and has diverse biological activities.¹⁻³ It is active against the bacteria *S. aureus*, *S. epidermidis*, *P. aeruginosa*, *E. cloacae*, *K. pneumoniae*, and *E. coli* (MICs = 0.6, 0.68, 0.97, 0.85, 1, and 1.08 mg/ml, respectively) and the fungi *C. albicans*, *C. tropicalis*, and *C. glabrata* (MICs = 1.85, 1.7, and 1.65 mg/ml, respectively).³ Androsin (10 mg/kg) inhibits bronchial obstruction in a guinea pig model of asthma induced by ovalbumin or platelet-activating factor (PAF).¹

References

1. Dorsch, W., Stuppner, H., Wagner, H., *et al.* Antiasthmatic effects of *Picrorhiza kurroa*: Androsin prevents allergen- and PAF-induced bronchial obstruction in guinea pigs. *Int. Arch. Allergy Appl. Immunol.* **95**(2-3), 128-133 (1991).
2. Stuppner, H. and Wagner, H. Minor iridoid and phenol glycosides of *Picrorhiza kurroa*. *Planta Med.* **55**(5), 467-469 (1988).
3. Zajdel, S.M., Graikou, K., Sotiroudis, G., *et al.* Two new iridoids from selected *Penstemon* species - antimicrobial activity. *Nat. Prod. Res.* **27**(24), 2263-2271 (2013).

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, 12/19/2022

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