

PRODUCT INFORMATION



Wilforgine

Item No. 35029

CAS Registry No.: 37239-47-7

Formal Name: 3-furancarboxylic acid, (8R,9R,10R,1S,12R,13R,14R,15S,18S,21S,22S,23R)-10,13,22,23-tetrakis(acetyloxy)-12-[(acetyloxy)methyl]-7,8,9,10,12,13,14,15,17,18,19,20-dodecahydro-21-hydroxy-8,18,21-trimethyl-5,17-dioxo-8,11-epoxy-9,12-ethano-11,15-methano-5H,11H-[1,9]dioxacyclooctadecino[4,3-b]pyridin-14-yl ester

MF: $C_{41}H_{47}NO_{19}$

FW: 857.8

Purity: $\geq 98\%$

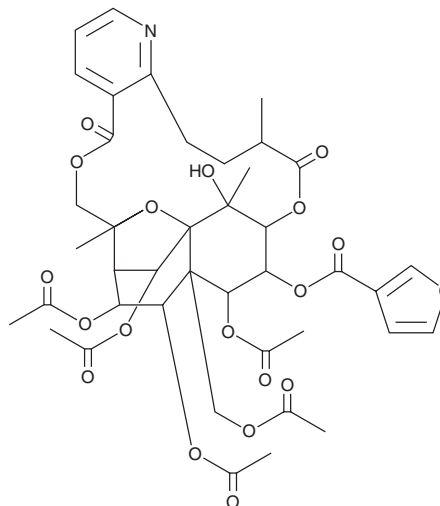
UV/Vis.: λ_{max} : 229 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Tripterygium wilfordii*



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

Laboratory Procedures

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

Description

Wilforgine is a sesquiterpene alkaloid that has been found in *T. wilfordii* and has paralytic activity.^{1,2} It induces paralysis in third instar *M. separata* larvae with a 50% narcotic concentration (NC_{50}) value of 0.0086 $\mu\text{mol/larvae}$.²

References

- Ouyang, X.-K., Jin, M.-C., and He, C.-H. Simultaneous determination of four sesquiterpene alkaloids in *Tripterygium wilfordii* Hook. F. extracts by high-performance liquid chromatography. *Phytochem. Anal.* **18**(4), 320-325 (2007).
- Qu, H., Lv, M., Yu, X., et al. Discovery of some piperine-based phenylsulfonylhydrazone derivatives as potent botanically narcotic agents. *Sci. Rep.* **5**, 13077 (2015).

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, 12/09/2022

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