

PRODUCT INFORMATION



Thermopsine

Item No. 33853

CAS Registry No.: 486-90-8
Formal Name: (7R,14R,14aS)-1,3,4,6,7,13,14,14a-octahydro-7,14-methano-2H,11H-dipyrido[1,2-a:1',2'-e][1,5]diazocin-11-one

Synonyms: NSC 21725, (-)-Thermopsine

MF: $C_{15}H_{20}N_2O$

FW: 244.3

Purity: $\geq 98\%$

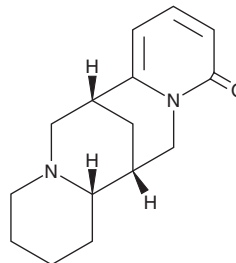
UV/Vis.: λ_{max} : 234, 311 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

Item Origin: Plant/*Paeonia lactiflora* Pall. root



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

Laboratory Procedures

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

Description

Thermopsine is an alkaloid that has been found in *Sophora tonkinensis* and has anti-inflammatory activity.¹ It reduces LPS-induced production of TNF- α in RAW 264.7 cells when used at a concentration of 50 μM .

Reference

1. He, L.-J., Liu, J.-S., Luo, D., *et al.* Quinolizidine alkaloids from *Sophora tonkinensis* and their anti-inflammatory activities. *Fitoterapia* **139**, 104391 (2019).

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, 10/25/2022

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