

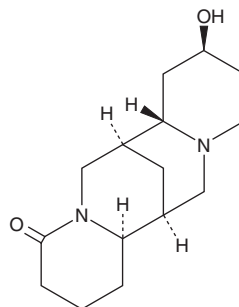
# PRODUCT INFORMATION



## Hydroxylupanine

Item No. 31171

**CAS Registry No.:** 15358-48-2  
**Formal Name:** (2S,7S,7aR,14S,14aS)-11H-dipyrido[1,2-a:1',2'-e][1,5]diazocin-11-one, dodecahydro-2-hydroxy-7,14-methano-2H  
**Synonyms:** (+)-Hydroxylupanine, 13 $\alpha$ -Hydroxylupanine  
**MF:** C<sub>15</sub>H<sub>24</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 264.4  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°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

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

### Description

Hydroxylupanine is an alkaloid that has been found in *L. lanatus*.<sup>1</sup> It is an inhibitor of ganglionic transmission.<sup>2</sup>

### References

1. Neto, A.T., Oliveira, C.Q., Ilha, V., *et al.* Quinolizidine alkaloids from *Lupinus lanatus*. *J. Mol. Struct.* **1004**(1-3), 174-177 (2011).
2. Petterson, D.S., Greirson, B.N., Allen, D.G., *et al.* Disposition of lupanine and 13-hydroxylupanine in man. *Xenobiotica* **24**(9), 933-941 (1994).

#### 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, 10/24/2022

#### 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](http://WWW.CAYMANCHEM.COM)