

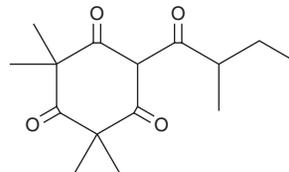
# PRODUCT INFORMATION



## Isoleptospermone

Item No. 44965

CAS Registry No.: 7375-66-8  
Formal Name: 2,2,4,4-tetramethyl-6-(2-methyl-1-oxobutyl)-1,3,5-cyclohexanetrione  
MF: C<sub>15</sub>H<sub>22</sub>O<sub>4</sub>  
FW: 266.3  
Purity: ≥95%  
Supplied as: An oil  
Storage: -20°C  
Stability: ≥2 years



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

### Laboratory Procedures

Isoleptospermone is supplied as an oil. A stock solution may be made by dissolving the isoleptospermone in the solvent of choice, which should be purged with an inert gas. Isoleptospermone is soluble in methanol and DMSO.

### Description

Isoleptospermone is a polyketide synthase-derived  $\beta$ -triketone that has been found in *L. scoparium* and has diverse biological activities.<sup>1-4</sup> It inhibits 4-hydroxyphenylpyruvate dioxygenase (HPPD; IC<sub>50</sub> = 14.3  $\mu$ M for the *A. thaliana* enzyme).<sup>1</sup> Isoleptospermone is active against *B. subtilis* and *T. mentagrophytes* (MICs = 2 and 1  $\mu$ g/ml, respectively).<sup>2</sup> It is lethal to *T. ni* and *C. includens* third-instar larvae (LC<sub>50</sub>s = 19.7 and 20.4  $\mu$ g/larva, respectively) and induces foliar damage in cabbage, soybean, and corn.<sup>3</sup> Isoleptospermone is also active against *S. scabiei* mites, young eggs, and mature eggs (LC<sub>50</sub>s = 59.4, 40.3, and 101.7 mM, respectively).<sup>4</sup>

### References

1. Dayan, F.E., Singh, N., McCurdy, C.R., *et al.*  $\beta$ -triketone inhibitors of plant p-hydroxyphenylpyruvate dioxygenase: Modeling and comparative molecular field analysis of their interactions. *J. Agric. Food Chem.* **57(12)**, 5194-5200 (2009).
2. van Klink, J.W., Larsen, L., Perry, N.B., *et al.* Triketones active against antibiotic-resistant Bacteria: synthesis, structure-activity relationships, and mode of action. *Bioorg. Med. Chem.* **13(24)**, 6651-6662 (2005).
3. Muturi, E.J., Neidmann, M.A., Ribeiro, V.P., *et al.* Insecticidal activity of manuka essential oil triketones against two major lepidopteran pests. *Pest Manag. Sci.* (2026).
4. Nammunige, N.A., Agnew-Francis, K.A., Fernando, D.D., *et al.*  $\beta$ -Triketones from *Leptospermum scoparium* (mānuka) oil show potential as scabicides. *Phytomedicine* **136**, 156321 (2025).

#### 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, 01/29/2026

#### 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