

PRODUCT INFORMATION



Bee Venom (*A. mellifera*)

Item No. 60500

Overview and Properties

Source: Bee venom (*Apis mellifera*)
Molecular Weight: 14 kDa
Storage: -20°C (as supplied)
Stability: ≥2 years
Supplied as: A lyophilized powder
Specific Activity: *batch specific* U/mg
Unit Definition: One unit of enzyme hydrolyzes one μmol of diheptanoyl thio-PC per minute at 25°C.
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

A. mellifera bee venom contains a variety of venom toxins, as well as proteins involved in antioxidant systems, protein folding, transport, carbohydrate metabolism, and energy metabolism.¹ Secretory phospholipase A₂ (sPLA₂) (Type III), also known as bee venom PLA₂ (bv-PLA₂), comprises 10-12% of crude bee venom dry weight and is the major bee venom toxin responsible for anaphylaxis.² However, bee venom and bv-PLA₂ also have anti-inflammatory and neuroprotective effects in rodent models.^{2,3}

References

1. Li, R., Zhang, L., Fang, Y., *et al.* Proteome and phosphoproteome analysis of honeybee (*Apis mellifera*) venom collected from electrical stimulation and manual extraction of the venom gland. *BMC Genomics* **14**, 766 (2013).
2. Lee, G. and Bae, H. Bee venom phospholipase A₂: Yesterday's enemy becomes today's friend. *Toxins (Basel)* **8(2)**, 48 (2016).
3. Zhang, S., Liu, Y., Ye, Y., *et al.* Bee venom therapy: Potential mechanisms and therapeutic applications. *Toxicon* **148**, 64-73 (2018).

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, 04/26/2024

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