

PRODUCT INFORMATION



Convulxin

Item No. 19082

CAS Registry No.: 37206-04-5

MW: ~84 kDa

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Hygroscopic, light sensitive, avoid freeze/thaw cycles

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

Description

Convulxin is a toxin that was first isolated from rattlesnake venom. It activates the collagen receptor glycoprotein VI (GPVI) on platelets and initiates blood clotting.^{1,2} Convulxin is used to study signaling through GPVI and to characterize platelet function.³⁻⁵

References

1. Vargaftig, B.B., Prado-Francheschi, J., Chignard, M., *et al.* Activation of guinea-pig platelets induced by convulxin, a substance extracted from the venom of *Crotalus durissus cascavella*. *Eur. J. Pharmacol.* **68(4)**, 451-464 (2016).
2. Jandrot-Perrus, M., Lagrue, A.-H., Okuma, M., *et al.* Adhesion and activation of human platelets induced by convulxin involve glycoprotein VI and integrin $\alpha 2\beta 1$. *J. Biol. Chem.* **272(43)**, 27035-27041 (1997).
3. Chatterjee, M.S., Denney, W.S., Jing, H., *et al.* Systems biology of coagulation initiation: Kinetics of thrombin generation in resting and activated human blood. *PLoS Comput. Biol.* **6(9)**, (2010).
4. Casari, C., Berrou, E., Lebret, M., *et al.* von Willebrand factor mutation promotes thrombocytopeny by inhibiting integrin $\alpha \text{IIb}\beta 3$. *J. Clin. Invest.* **123(12)**, 5071-5081 (2013).
5. van Bladel, E.R., Laarhoven, A.G., van der Heijden, L.B., *et al.* Functional platelet defects in children with severe chronic ITP as tested with 2 novel assays applicable for low platelet counts. *Blood* **123(10)**, 1556-1563 (2014).

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, 12/10/2025

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