

# Charybdotoxin, (recombinant)

**K<sup>+</sup> channel blocker**

Charybdotoxin was originally isolated from the venom of the Israeli scorpion *Leiurus quinquestriatus hebraeus*. Charybdotoxin blocks K<sub>Ca</sub>1.1 (large conductance Ca<sup>2+</sup> activated K<sup>+</sup>, Slo) channels in nanomolar concentrations as well as K<sub>V</sub>1.2 (K<sub>d</sub>, 14 nM) and K<sub>V</sub>1.3 (K<sub>d</sub>, 2.6 nM) channels.). However, experiments with cloned K<sub>Ca</sub>1.1 channels demonstrate the strong effect of the slo subunits on the potency of block by Charybdotoxin. This product is a recombinant peptide expressed in and extracted from *E. coli*.

## Ordering Information

[Order Online »](#)

BML-KC168-0100	100µg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Appearance      Lyophilized solid.

CAS      95751-30-7

MW      4313

Purity      ≥98%

Sequence      Gln-Phe-Thr-Asn-Val-Ser-Cys-Thr-Thr-Ser-Lys-Glu-Cys-Trp-Ser-Val-Cys-Gln-Arg-Leu-His-Asn-Thr-Ser-Arg-Gly-Lys-Cys-Met-Asn-Lys-Lys-Cys-Arg-Cys-Tyr-Ser

Solubility      Soluble in Water.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)