

PRODUCT INFORMATION



Myr-Tat-CBD3 (trifluoroacetate salt)

Item No. 37465

Formal Name:	N-myristoyl-L-tyrosylglycyl-L-arginyl-L-lysyl-L-lysyl-L-arginyl-L-arginyl-L-glutamyl-L-arginyl-L-arginyl-L-arginyl-L-alanyl-L-arginyl-L-seryl-L-arginyl-L-leucyl-L-alanyl-L- α -glutamyl-L-leucyl-L-arginylglycyl-L-valyl-L-prolyl-L-arginylglycyl-L-leucine, trifluoroacetate salt	Myr – Tyr – Gly – Arg – Lys – Lys – Arg – Arg – Gln – Arg – Arg – Arg – Ala – Arg – Ser – Arg – Leu – Ala – Glu – Leu – Arg – Gly – Val – Pro – Arg – Gly – Leu – OH
Synonyms:	N-myristate-Tat-CBD3, Myr-Tat-Calcium Channel-binding Domain 3	
Peptide Sequence:	myristoyl-YGRKKRRQRRRARSRLAELRGVPRGL-OH	
MF:	$C_{148}H_{269}N_{59}O_{33} \cdot XCF_3COOH$	$\bullet XCF_3COOH$
FW:	3,403.1	
Purity:	$\geq 95\%$	
Supplied as:	A solid	
Storage:	-20°C	
Stability:	≥ 4 years	

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

Laboratory Procedures

Myr-Tat-CBD3 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Myr-Tat-CBD3 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Myr-Tat-CBD3 is an inhibitor of the protein-protein interaction between the N-type voltage-gated calcium channel $Ca_v2.2$ and collapsin response mediator protein 2 (CRMP2).¹ It inhibits the $Ca_v2.2$ -CRMP2 interaction by approximately 81% when used at a concentration of 10 μM and inhibits potassium-induced calcium influx in primary rat dorsal root ganglion (DRG) neurons ($IC_{50} = \sim 2.8 \mu M$). Myr-Tat-CBD3 (20 μM) inhibits calcium, but not sodium, currents in primary DRG neurons. Intrathecal administration of myr-Tat-CBD3 (20 $\mu g/5 \mu l$) prevents carrageenan-induced decreases in paw withdrawal latencies in rats. Myr-Tat-CBD3 decreases cue-induced reinstatement of cocaine-seeking behavior in rats.²

References

1. François-Moutal, L., Wang, Y., Moutal, A., *et al.* A membrane-delimited N-myristoylated CRMP2 peptide aptamer inhibits $Ca_v2.2$ trafficking and reverses inflammatory and postoperative pain behaviors. *Pain* **156(7)**, 1247-1264 (2015).
2. Buchta, W.C., Moutal, A., Hines, B., *et al.* Dynamic CRMP2 regulation of $Ca_v2.2$ in the prefrontal cortex contributes to the reinstatement of cocaine seeking. *Mol. Neurobiol.* **57(1)**, 346-357 (2020).

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, 10/09/2023

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