

PRODUCT INFORMATION



Tat-AKAP79 (326-336) (trifluoroacetate salt)

Item No. 37444

| | | |
|--------------------------|---|--|
| Synonym: | Tat-A-kinase Anchor Protein 79 (326-336) | |
| Peptide Sequence: | YGRKKRRQRRRKSKSEESKRME-OH | H-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg- |
| MF: | $C_{117}H_{211}N_{49}O_{34}S \cdot XCF_3COOH$ | Arg-Lys-Ser-Lys-Ser-Glu-Glu-Ser-Lys-Arg- |
| FW: | 2,880.3 | Met-Glu-OH |
| Purity: | ≥98% | • XCF ₃ COOH |
| 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

Tat-AKAP79 (326-336) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Tat-AKAP79 (326-336) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Tat-AKAP79 (326-336) is a peptide composed of the HIV-1 Tat protein transduction domain linked to an 11-amino acid peptide corresponding to residues 326-336 of A-kinase anchor protein 79 (AKAP79).¹ It inhibits protein kinase A (PKA) or PKC activation-induced sensitization of transient receptor potential vanilloid 1 (TRPV1) in isolated rat dorsal root ganglion (DRG) neurons when used at a concentration of 200 μM. *In vivo*, Tat-AKAP79 (326-336) inhibits nocifensive behavior induced by formalin or phorbol 12-myristate 13-acetate (PMA; Item No. 10008014) in mice. It also increases the latency to paw withdrawal from a radiant heat stimulus and increases the mechanical threshold force to paw withdrawal in a mouse model of carrageenan-induced inflammatory pain.

Reference

1. Btsh, J., Fischer, M.J.M., Stott, K., *et al.* Mapping the binding site of TRPV1 on AKAP79: Implications for inflammatory hyperalgesia. *J. Neurosci.* **33(21)**, 9184-9193 (2013).

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, 10/11/2023

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