

PRODUCT INFORMATION



Tat (48-60) (trifluoroacetate salt)

Item No. 37499

Formal Name: glycyl-L-arginyl-L-lysyl-L-lysyl-L-arginyl-L-arginyl-L-glutamyl-L-arginyl-L-arginyl-L-arginyl-L-prolyl-L-prolyl-L-glutamine, trifluoroacetate salt

Synonym: Transactivator of Transcription Peptide (48-60) H—Gly—Arg—Lys—Lys—Arg—Arg—Gln—Arg—Arg—Arg—Pro—Pro—Gln—OH

Peptide Sequence: H-GRKKRRQRRRPPQ-OH

MF: $C_{70}H_{131}N_{35}O_{16} \cdot XCF_3COOH$

FW: 1,719.0

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

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

Laboratory Procedures

Tat (48-60) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Tat (48-60) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Tat (48-60) (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Tat (48-60) (trifluoroacetate salt) in these solvents is approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Tat (48-60) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Tat (48-60) (trifluoroacetate salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Tat (48-60) is an arginine-rich peptide, a fragment of HIV Tat, and a form of the cell-penetrating peptide Tat (48-60) amide.

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, 10/05/2023

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