

# PRODUCT INFORMATION



## Tat (2-4) (trifluoroacetate salt)

Item No. 37445

**Synonym:** Transactivator of Transcription Peptide (2-4)  
**Peptide Sequence:** YGRKKRRQRRRGYGRKKRRQRRRG-OH      H-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-  
**MF:**  $C_{132}H_{240}N_{66}O_{29} \cdot XCF_3COOH$       Arg-Gly-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-  
**FW:** 3,215.8      Arg-Arg-Arg-Gly-OH  
**Purity:**  $\geq 98\%$       •  $XCF_3COOH$   
**Supplied as:** A solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years

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

### Laboratory Procedures

Tat (2-4) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Tat (2-4) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Tat (2-4) is a peptide derived from HIV-1 transactivator of transcription (Tat) protein. It is composed of two repeats of the arginine-rich cationic peptide Tat (47-58).

#### 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, 10/04/2023

#### 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