

# PRODUCT INFORMATION



## Tat (47-57)

Item No. 43817

CAS Registry No.: 191936-91-1

Formal Name: L-tyrosylglycyl-L-arginyl-L-lysyl-L-lysyl-L-arginyl-L-arginyl-L-glutamyl-L-arginyl-L-arginyl-L-arginine

H-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-

Peptide Sequence: YGRKKRRQRRR-OH

Arg-OH

Synonyms: Tat11,

Transactivator of Transcription Peptide (47-57)

MF:  $C_{64}H_{118}N_{32}O_{14}$

FW: 1,559.8

Purity:  $\geq 95\%$

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 (47-57) is supplied as a solid. A stock solution may be made by dissolving the Tat (47-57) in the solvent of choice, which should be purged with an inert gas. Tat (47-57) is slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

Tat (47-57) is an arginine-rich cell-penetrating peptide (CPP) and a synthetic fragment of HIV-1 Tat.<sup>1,2</sup> It has been used to deliver different types of cargo to cells, including fluorescent proteins and small molecules.

### References

1. Patel, S.G., Sayers, E.J., He, L., *et al.* Cell-penetrating peptide sequence and modification dependent uptake and subcellular distribution of green fluorescent protein in different cell lines. *Sci. Rep.* **9(1)**, 6298 (2019).
2. Goswami, D., Vitorino, H.A., Alta, R.Y., *et al.* Deferasirox-TAT(47-57) peptide conjugate as a water soluble, bifunctional iron chelator with potential use in neuromedicine. *Biometals* **28(5)**, 869-877 (2015).

#### 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, 10/07/2025

#### 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