



## Product Data Sheet

---

<b>Product Name:</b>	TAT-NSF700 Fusion Peptide	
<b>Catalog Number:</b>	AS-62238 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-Arg-Gly-Gly-Gly-Leu-Leu-Asp-Tyr-Val-Pro-Ile-Gly-Pro-Arg-Phe-Ser-Asn-Leu-Val-Leu-Gln-Ala-Leu-Leu-Val-Leu-OH (3-letter code) YGRKKRRQRRR-GGG-LLDYVPIGPRFSNLVLQALLVL (1-letter code)	
<b>Molecular Weight:</b>	4167.0	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** TAT-NSF700 Fusion peptide is freely soluble in water.

**Storage:** TAT-NSF700 Fusion peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** This peptide is the N-Ethyl-maleimide-sensitive factor (NSF) inhibitor fusion polypeptide composed of 11-amino acid cell permeable HIV transactivating regulatory protein (TAT) domain fused to a 22 amino acid NSF domain. TAT-NSF700 inhibits thrombin-induced exocytosis of endothelial cells in a dose-responsive manner. Ref: Matsushita, K. et al. *Mol. Pharmacol.* **67**,1137 (2005).

**Additional Information:** Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

The control group was treated with a fusion peptide possessing identical amino acid composition but with a scrambled sequence TAT-NSF700scr (5  $\mu\text{M}$ ) 60 min before ATP stimulation. TAT-NSF fusion peptides were obtained from Anaspec (1000x PBS)- [Trang, T. et al. \*J Neurosci\* \*\*29\*\*, 3518 \(2009\).](#)

### Published Citations:

Matsushita, K. et al. *Mol Pharmacol* **67**, 1137 (2005).  
Trang, T. et al. *J Neurosci* **29**, 3518 (2009).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
TAT-NSF222scr Fusion Polypeptide, scrambled	AS-62210	1 mg
YGRKKRRQRRR-GGG-ENSFRFLADIFPAKAFPVRFE		

TAT-NSF222 Fusion Peptide YGRKKRRQRRR-GGG-LDKEFNSIFRRAFASRVFPPE	AS-62240	1 mg
Cys(Npys)-TAT (47-57), FAM-labeled C(Npys)YGRKKRRQRRR-K(FAM)-NH <sub>2</sub>	AS-61213	1 mg

*For Research Use Only*