

Product Data Sheet

Product Name: Cys-TAT (47-57)

Catalog Number: AS-61212 (1 mg) Lot Number: See label on vial

Sequence: H-Cys-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-Arg-NH2 (3-letter code)

CYGRKKRRQRRR-NH2 (1-letter code)

Molecular Weight: 1662.0

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Cys-TAT (47-57) peptide is freely soluble in water.

Storage: Cys-TAT (47-57) peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20 °C or lower.

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.

Peptides were synthesized and conjugated to TAT carrier peptide (residues 47-57) via cysteine S-S bond by Anaspec, San Jose, CA. Hearts were submerged into a heat-jacketed organ bath at 37 °C. Coronary effluent was collected to determine creatine phosphokinase release. After 10 min of equilibration, the hearts were subjected to 40 min of global ischemia and 60 min of reperfusion. The hearts were perfused with 1 µm TAT-conjugated peptide for 10 min prior to ischemia. In addition to the relative amount of creatine phosphokinase released, a measure of cardiac myocyte lysis, tissue damage was assessed by triphenyltetrazolium chloride (TTC) staining of heart crosssections to quantitate the amount of infarcted (dead) tissue-Brandman, R. et al. J. Biol. Chem. 282, 4113 (2007).

Published Citations:

Brandman, R. et al. J. Biol. Chem. 282, 4113 (2007).

Related Products:

Name	Cat #	Size
TAT (47-57)	AS-60023-1	1 mg
YGRKKRRQRRR	AS-60023-5	5 mg
TAT (47-57), FAM-labeled FAM-YGRKKRRQRRR	AS-61210	1 mg

TAT (47-57), TAMRA-labeled TAMRA-YGRKKRRQRRR	AS-61211	1 mg
FITC-LC-TAT (47-57) FITC-LC-YGRKKRRQRRR-NH2	AS-27042 AS-27043	0.5 mg 1 mg
TAT (47-57) GGG-Cys(Npys) YGRKKRRQRRRGGG-C(Npys)-NH2	AS-61215	1 mg
Biotin-TAT (47-57) Biotin-YGRKKRRQRRR	AS-61209	1 mg
Cys(Npys)-TAT (47-57), FAM-labeled C(Npys)YGRKKRRQRRR-K(FAM)-NH2	AS-61213	1 mg
Cys(Npys)-TAT (47-57) C(Npys)YGRKKRRQRRR-NH2	AS-61214	1 mg

For Research Use Only