



Product Data Sheet

Product Name: T20
Catalog Number: 61235 (1 mg) **Lot Number:** See label on vial
Sequence: Ac-Tyr-Thr-Ser-Leu-Ile-His-Ser-Leu-Ile-Glu-Glu-Ser-Gln-Asn-Gln-Gln-Glu-Lys-Asn-Glu-Gln-Glu-Leu-Leu-Glu-Leu-Asp-Lys-Trp-Ala-Ser-Leu-Trp-Asn-Trp-Phe-NH₂ (3-letter code)
Ac-YTSLIHSLIEESQNQQEKNEQELLELDKWASLWNWF-NH₂ (1-letter code)
Molecular Weight: 4493.0
% Peak Area by HPLC: ≥ 95
Appearance: Lyophilized white powder

Peptide Reconstitution: Reconstitute by adding 100 µl 1% NH₄OH to 1 mg T20 peptide. Dilute this peptide solution to approximately 1 mg/ml (or more dilute) with a buffer such as PBS or another buffer; aliquot and store at -20°C.

Storage: T20 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.

Description: This 36-residue peptide corresponds to aa 643-678 of the C-terminus of HIV-1LAIgp41. It strongly inhibits HIV-1 viral fusion with an EC₅₀ of 1 ng/ml. Ref: Lawless, MK. et al. *Biochem.* **35**, 13697 (1996); Su, SB. et al. *Blood* **93**, 3885 (1999).

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 cells were exposed to serial three-fold dilutions of cell-free supernatants harvested from the cotransfected cells containing 40 µg/ml DEAE-dextran (Sigma) in 96-well plates. After 6 h exposure at 37 °C in an atmosphere of 5% CO₂, the viral fusion inhibitor T-20 (AnaSpec, San Jose, CA) was added at a final concentration of 2 µg/ml. After 48 h of incubation, 100 µl of the Bright-Glo luciferase assay system (Promega) was added to each well. The plates were read in a Wallac Victor2 1420 multilabel counter (Perkin-Elmer) for luciferase activity. All viral infection assays were performed in triplicate-[Li, PL. et al. *Virology* **331**, 367 \(2005\).](#)

Published Citations:

Li, PL. et al. *Virology* **331**, 367 (2005).

Related Products:

Name	Cat #	Size
T20-Lys(FAM) (Ac-YTSLIHSLIEESQNQQEKNEQELLELDKWASLWNWF-K(FAM)-NH ₂)	61236	0.1 mg
T20-Lys(TAMRA) (Ac-YTSLIHSLIEESQNQQEKNEQELLELDKWASLWNWF-K(TAMRA)-NH ₂)	61237	0.1 mg
T20-Lys(HiLyte Fluor™488) (Ac-YTSLIHSLIEESQNQQEKNEQELLELDKWASLWNWF-K(HiLyte Fluor™488)-NH ₂)	61238	0.1 mg

For Research Use Only