

HIV-1 TAT Clade-C

Ordering Information

[Order Online »](#)

| | |
|-------------|------|
| JBS-PR-1206 | 10µg |
|-------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Human Immunodeficiency Virus 1 Antigen

Formulation lyophilised (with no additives)

MW 21 kDa

Purity > 90 % (SDS-PAGE, HPLC)

Solubility It is recommended to reconstitute the lyophilised HIV-1 TAT in bidest H₂O not less than 100 µg/ml, which can then be further diluted to other aqueous solutions. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).

Source recombinant, E. coli

Specificity Reacts with anti-Tat polyclonal antibodies from human, monkey, rabbit and mouse serum. Recognized by anti-Tat (HIV-1) polyclonal antibody.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com