

ab218268 HIV-1 p24 SimpleStep ELISA® Kit

For the quantitative measurement of HIV-1 p24 protein in human serum, plasma, cell culture supernatant, and cell and tissue extract samples.

For overview, typical data and additional information please visit: www.abcam.com/ab218268

Storage and Stability: Store kit at +4°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted. Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

Materials Supplied

Item	Quantity	Storage Condition
HIV-1 p24 Capture Antibody 10X	600 µL	+4°C
HIV-1 p24 Detector Antibody 10X	600 µL	+4°C
HIV-1 p24 Lyophilized Recombinant Protein	2 Vials	+4°C
Antibody Diluent 5BI	6 mL	+4°C
10X Wash Buffer PT	20 mL	+4°C
Cell Extraction Buffer PTR 5X	10 mL	+4°C
TMB Substrate	12 mL	+4°C
Stop Solution	12 mL	+4°C
Sample Diluent NS	50 mL	+4°C
Sample Diluent 50BS	20 mL	+4°C
Anti-tag coated microplate (12 x 8 well strips)	96 wells	+4°C
Plate Seal	1	+4°C
50X Cell Extraction Enhancer Solution *	1 mL	+4°C

***50X Cell Extraction Enhancer Solution is included in the kit but is not used for testing any currently validated sample types**

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

Microplate reader capable of measuring absorbance at 450 or 600 nm.
 Method for determining protein concentration (BCA assay recommended).
 Deionized water.
 Multi- and single-channel pipettes.
 Tubes for standard dilution.
 Plate shaker for all incubation steps.
 Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).
 1N HCL, 1.2 N NaOH.

Reagent Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells. The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.

Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations

1X Cell Extraction Buffer PTR (For cell and tissue extracts only) Prepare 1X Cell Extraction Buffer PTR by diluting Cell Extraction Buffer PTR 5X to 1X with deionized water. To make 10 mL 1X Cell

Extraction Buffer PTR combine 8 mL deionized water and 2 mL Cell Extraction Buffer PTR 5X. Mix thoroughly and gently. If required protease inhibitors can be added.

1X Wash Buffer PT Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL Wash Buffer PT 10X with 45 mL deionized water. Mix thoroughly and gently.

Antibody Cocktail: Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 5BI. To make 3 mL of the Antibody Cocktail combine 300 µL 10X Capture Antibody and 300 µL 10X Detector Antibody with 2.4 mL Antibody Diluent 5BI. Mix thoroughly and gently.

Standard Preparation

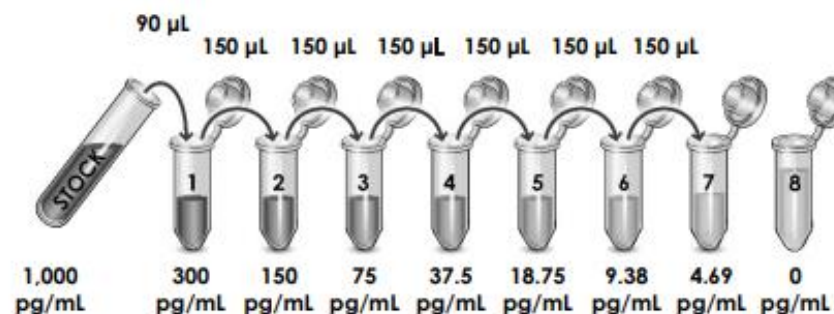
Always prepare a fresh set of standards for every use. Discard working standard dilutions after use.

The following section describes the preparation of a standard curve for duplicate measurements (recommended).

The standard of this kit corresponds in amino acids sequence to HIV-1 p24, group M, subtype B, strain 92418.

IMPORTANT: If the protein standard vial has a volume identified on the label, reconstitute the HIV-1 p24 standard by adding that volume of Diluent indicated on the label. Alternatively, if the vial has a mass identified, reconstitute the HIV-1 p24 standard by adding 500 µL Diluent. Hold at room temperature for 10 minutes and mix gently. This is the 1,000 pg/mL **Stock Standard** Solution.

- For serum and plasma samples measurements**, reconstitute the HIV-1 p24 standard by adding Sample Diluent 50BS.
For cell culture supernatant samples reconstitute the HIV-1 p24 standard by adding Sample Diluent NS
For cell and tissue extract samples reconstitute the HIV-1 p24 standard by adding 1X Cell Extraction Buffer PTR.
- Label eight tubes, Standards 1–8.
- Add 210 µL of appropriate diluent to tube number 1 and 150 µL of appropriate diluent into numbers 2-8.
- Use the Stock Standard to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:



To convert sample values obtained with the kit to approximate NIBSC 90/636 units, use the following equation: NIBSC 90/636 approximate value (IU/mL) = 0.43 x SimpleStep HIV-1 p24 value (pg/mL).

Sample Preparation

Plasma: Collect plasma using EDTA or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Samples may be assayed neat or dilute samples into Sample Diluent 50BS and assay. Store undiluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles. Acid Treatment Protocol: For the dissociation of immune complex in plasma samples, add 150 µL of sample + 75 µL of 1N HCL and mix. Incubate 10 minutes at room temperature. Add 75 µL of 1.2 N NaOH + 0.5 M HEPES and mix. Final sample dilutions are 1:2.

Serum: Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 2,000 x g for 10 minutes and collect serum. Samples may be assayed neat or dilute samples into Sample Diluent 50BS and assay. Store undiluted serum at -20°C or below. Avoid repeated freeze-thaw cycles. Acid Treatment Protocol: For the dissociation of immune complex in serum samples, add 150 µL of sample + 75 µL of 1N HCL and mix. Incubate 10 minutes at room temperature. Add 75 µL of 1.2 N NaOH + 0.5 M HEPES and mix. Final sample dilutions are 1:2.

Cell Culture Supernatants: Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants and assay or dilute samples into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. Avoid repeated freeze-thaw cycles.

Preparation of extracts from cell pellets: Collect non-adherent cells by centrifugation or scrape to collect adherent cells from the culture flask. Typical centrifugation conditions for cells are 500 x g for 5 minutes at 4°C. Rinse cells twice with PBS. Solubilize pellet at 2x10⁷ cell/mL in chilled 1X Cell Extraction Buffer PTR. Incubate on ice for 20 minutes. Centrifuge at 18,000 x g for 20 minutes at 4°C. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay. Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

Preparation of extracts from adherent cells by direct lysis (alternative protocol): Remove growth media and rinse adherent cells 2 times in PBS. Solubilize the cells by addition of chilled 1X Cell Extraction Buffer PTR directly to the plate (use 750 µL - 1.5 mL 1X Cell Extraction Buffer PTR per confluent 15 cm diameter plate). Scrape the cells into a microfuge tube and incubate the lysate on ice for 15 minutes. Centrifuge at 18,000 x g for 20 minutes at 4°C. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay. Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

Preparation of extracts from tissue homogenates: Tissue lysates are typically prepared by homogenization of tissue that is first minced and thoroughly rinsed in PBS to remove blood (dounce homogenizer recommended). Homogenize 100 to 200 mg of wet tissue in 500 µL - 1 mL of chilled 1X Cell Extraction Buffer PTR. For lower amounts of tissue adjust volumes accordingly. Incubate on ice for 20 minutes. Centrifuge at 18,000 x g for 20 minutes at 4°C. Transfer the supernatants into clean tubes and discard the pellets. Assay samples immediately or aliquot and store at -80°C. The sample protein concentration in the extract may be quantified using a protein assay. Dilute samples to desired concentration in 1X Cell Extraction Buffer PTR.

Plate Preparation

The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents. Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.

For each assay performed, a minimum of two wells must be used as the zero control. For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates). Differences in well absorbance or "edge effects" have not been observed with this assay.

Assay Procedure

Equilibrate all materials and prepared reagents to room temperature prior to use. We recommend that you assay all standards, controls and samples in duplicate.

- 1) Prepare all reagents, working standards, and samples as directed in the previous sections.
- 2) Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, resealed and return to 4°C storage.
- 3) Add 50 µL of all sample or standard to appropriate wells.
- 4) Add 50 µL of the Antibody Cocktail to each well.
- 5) Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
- 6) Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and blot it against clean paper towels to remove excess liquid.
- 7) Add 100 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm. *Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes. Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.*
- 8) Add 100 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.
- 9) Alternative to 7 – 8: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

Mode	Kinetic
Wavelength	600 nm
Time	up to 20 min
Interval	20 sec - 1 min
Shake	Shake between readings

Note that an endpoint reading can also be recorded at the completion of the kinetic read by adding 100 µL Stop Solution to each well and recording the OD at 450 nm.

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

www.abcam.com/protocols/the-complete-elisa-guide

For technical support contact information, visit: www.abcam.com/contactus