

# Lentivirus Titer p24 ELISA Detection Kit

## Overview

Cat.No. HG-P001L

This product uses a double antibody sandwich method to detect HIV 1 p24 protein in samples. A monoclonal antibody specific to HIV 1 p24 antigen is coated on a microplate, and the standard or test sample is added into the reaction well. At the same time, the anti HIV 1 p24 secondary antibody is added and incubated at room temperature to form the antibody antigen secondary antibody complex. The unconjugated compounds are removed by washing and protein content in the sample is indicated by the intensity of TMB color development.

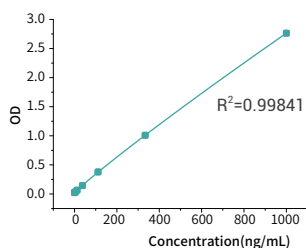
## Specification

**Assay range:** 1.37-1000 ng/mL

**Detection sensitivity:** 0.35 ng/mL

**Precision:** CV%≤10%, RE%≤±15%

## Standard curve



## Datasheet

Concentration of standard (ng/mL)	OD value1	OD value2	Mean value
200	2.846	2.841	2.844
100	1.965	1.979	1.972
50	1.110	1.189	1.150
25	0.636	0.582	0.609
12.5	0.362	0.338	0.350
6.25	0.213	0.204	0.209
0	0.061	0.059	0.060
0	0.022	0.022	0.022