

FOR RESEARCH USE ONLY!

Instant Lentivirus Detection Card

05/20

(Catalog # K1470-10, -20; 10 tests, 20 tests; Store at RT)

I. Introduction:

BioVision's **Instant Lentivirus Detection Card** is a rapid detection card which targets lentiviral capsule proteins as a marker to determine viral titer within 10 min. It is a simple lateral flow immunoassay and the results obtained are comparable to conventional infection-based assays that would take 3-4 days to complete. Results are in a red colour-readout format. The appearance of a control red band indicates a positive assay result. The card is recommended for detecting useable viral titer (at least 1×10^6 IU/ml) for downstream applications and/or to determine when to harvest lentiviral supernatant for optimal yield.

II. Application:

- A rapid tool to detect Lentivirus

III. Key Features:

- Rapid, easy and convenient
- Obtain results in 10-15 min
- Direct observation without any instruments
- **High Sensitivity**

IV. Sample Type:

Lentivirus packaging supernatant

V. Contents:

Components	K1470-10 (10 Tests)	K1470-20 (20 Tests)	Part Number
Detection Card	10	20	K1470-XX-1

VI. Shipping and Storage Conditions:

The card should be stored at room temperature upon arrival in the original packaging. Do not freeze the detection cards. The product is stable for up to 12 months when stored properly.

VII. Detection Protocol:

1. Place the **Detection card** horizontally on a level surface.
2. Add 100 μ l of **packaging supernatant** directly to the loading well.
3. Wait for 10-15 min until **red bands** appear. Observe the appearance of a red control band near the top of the card to determine a successful run. The intensity of the red band correlates with the relative viral titer.

Note: If no red control band shows up, the assay result is void.

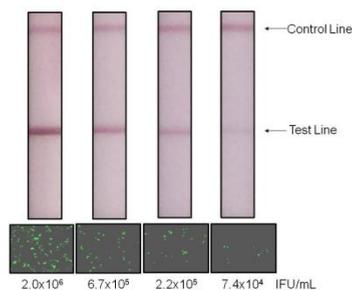


Figure shows the appearance of red bands for different dilutions of Lentivirus using Instant Lentivirus Detection Card

IX. Related Products:

BioVision Product Name	Cat. No.	Sizes
Mag-Lentivirus and Retrovirus Purification Kit	K1458	20, 100 Preps
Lentivirus Mini Purification Kit	K1305	-10, -20 preps
Lentivirus Maxi Purification Kit	K1306	-2, -4, -10 preps
PEG Virus Precipitation Kit	K904	-50, -200 preps
Mag-Adenovirus Purification Kit	K1459	-10, -200 preps
Adenovirus Mini Purification Kit	K1300	-10, -20 preps
Adenovirus Maxi Purification Kit	K1301	-2, -4, -10 preps
Retrovirus Mini Purification Kit	K1307	-10, -20 preps
Retrovirus Maxi Purification Kit	K1308	-2, -4, -10 preps

FOR RESEARCH USE ONLY! Not to be used on humans.