

# RCL(VSVG) Gene Copy Number Detection Kit (qPCR)

## Overview

Cat.No. HG-RC001

This kit is designed for the quantitative detection of RCL gene copy number in the genome of CAR-T cells prepared by using HIV-1 lentiviral vector technology.

This kit adopts the fluorescent probe method and multiplex PCR method to detect the DNA sequence related to integration or expression function on the transfer plasmid, and the VSVG gene copy number in the sample can be calculated. The kit is a rapid, specific and reliable device.

## Specification

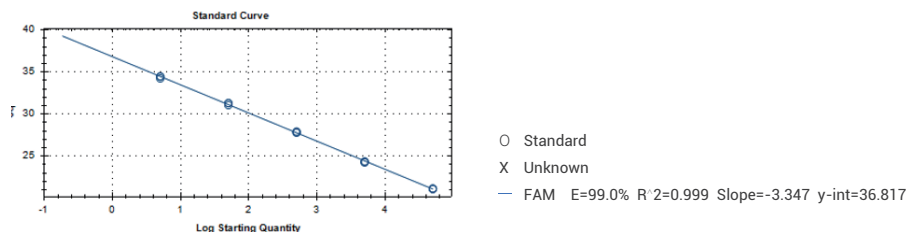
**Assay range:**  $1.00 \times 10^1 \sim 1.00 \times 10^5$  copies/ $\mu$ L

**Limit of quantitation:** 8 copies/ $\mu$ L

**Limit of detection:** 1 copies/ $\mu$ L

**Precision:** CV% $\leq$ 15%

## Standard curve



## Datasheet

Concentration (copies/ $\mu$ L)	Log10 Concentration	Ct Value(1)	Ct Value(2)	Ct Value(3)	Ct Mean Value	CV
1.00E+05	5.00	20.27	20.16	20.13	20.18	0.36%
1.00E+04	4.00	23.35	23.21	23.27	23.28	0.32%
1.00E+03	3.00	26.93	26.85	26.96	26.91	0.21%
1.00E+02	2.00	30.31	30.31	30.20	30.27	0.21%
1.00E+01	1.00	33.54	33.30	33.19	33.34	0.53%
Amplification efficiency						99.60%