

Vero Residual DNA Detection Kit (qPCR)

Overview

Cat.No. HG-VE001

This kit is designed for the quantitative detection of residual Vero host cell DNA in intermediates, semi-finished products and finished products of various biological products.

This kit adopts the principle of Taqman probe to quantitatively detect residual Vero DNA in samples. The kit is a rapid, specific and reliable device, with the minimum detection limit reaching fg level.

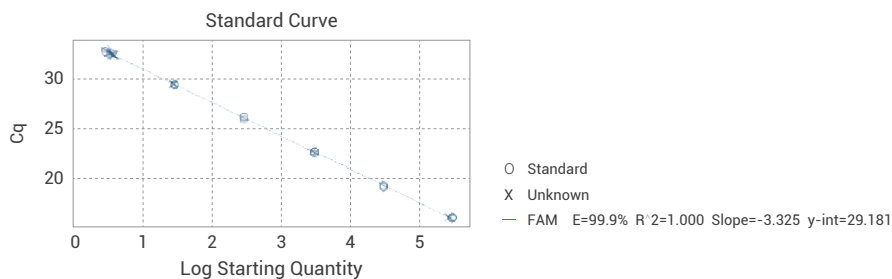
Specification

Assay range: 3.00~3.00x10⁵fg/μL

Limit of Quantitation: 3fg/μL

Precision: CV%≤15%

Standard curve



Datasheet

Concentration (fg/μl)	Log10 Concentration	Ct Value(1)	Ct Value(2)	Ct Value(3)	Ct Mean Value	CV
3.00E+05	5.48	11.07	10.99	11.15	11.07	0.72%
3.00E+04	4.48	14.09	14.18	14.31	14.20	0.78%
3.00E+03	3.48	17.68	17.46	17.64	17.59	0.65%
3.00E+02	2.48	20.76	21.07	21.03	20.95	0.80%
3.00E+01	1.48	24.32	24.19	24.22	24.24	0.29%
3.00E+00	0.48	27.84	27.55	27.54	27.64	0.61%
Amplification efficiency						99.9%