

# Human Residual DNA Detection Kit (qPCR)

## Overview

Cat.No. HG-HD001

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

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

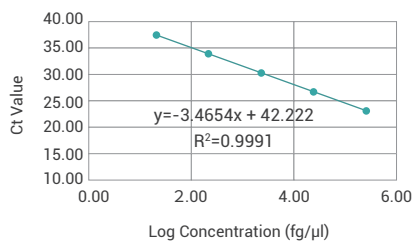
## Specification

**Assay range:**  $3.00 \times 10^1 \sim 3.00 \times 10^5$  fg/ $\mu$ L

**Limit of quantitation:**  $3.00 \times 10^1$  fg/ $\mu$ L

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

## Standard curve



## Datasheet

Concentration (fg/ $\mu$ l)	Log Concentration	Ct Value(1)	Ct Value(2)	Ct Value(3)	Ct Mean Value	Recovery rate
3.00E+05	5.48	23.52	23.37	23.39	23.43	88%
3.00E+04	4.48	26.64	26.55	26.45	26.54	111%
3.00E+03	3.48	30.16	30.07	30.06	30.09	105%
3.00E+02	2.48	33.75	33.33	33.52	33.53	107%
3.00E+01	1.48	38.3	36.49	36.90	37.26	90%
Amplification efficiency						94.34%