

E.coli Residual DNA Detection Kit (qPCR)

Overview

Cat.No. HG-ED001

This kit is designed for the quantitative detection of E.coli 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 E.coli residual DNA in samples. The kit is a rapid, specific and reliable device, with the minimum detection limit reaching fg level.

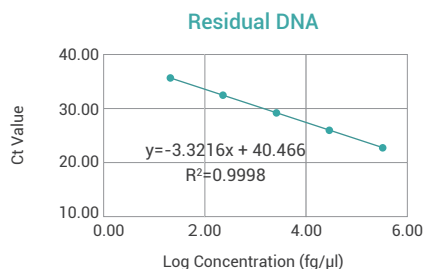
Specification

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

Limit of quantitation: 3.00×10^1 fg/ μ L

Precision: CV% \leq 15%

Standard curve



Datasheet

Concentration (fg/ μ l)	Log Concentration	Ct Value(1)	Ct Value(2)	Ct Value(3)	Ct Mean Value	Recovery rate
3.00E+05	5.48	22.30	22.33	22.32	22.32	97%
3.00E+04	4.48	25.51	25.43	25.49	25.48	108%
3.00E+03	3.48	29.09	28.89	28.90	28.96	97%
3.00E+02	2.48	32.38	32.30	32.72	32.32	95%
3.00E+01	1.48	35.52	35.28	35.72	35.51	104%
Amplification efficiency						100.01%