

E.coli Residual Total RNA Detection Kit (RT-PCR)

Overview

Cat.No. HG-ER001

This kit is designed for the quantitative detection of residual E.coli total RNA in various biological products to improve control quality of nucleic acid.

This kit adopts the principle of the RT-PCR fluorescent probe, combining reverse transcription PCR technology and fluorescent probe method, to realize one-step quantitative detection.

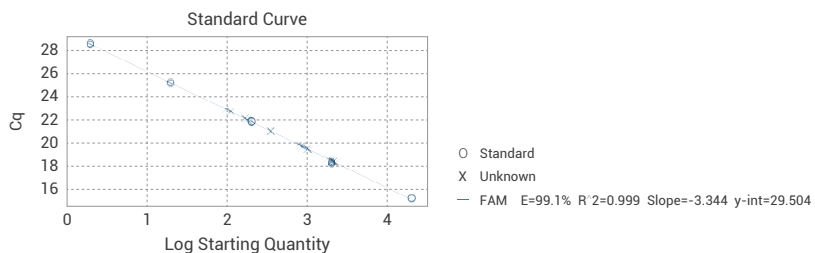
Specification

Assay range: $2.00 \times 10^{-3} \sim 2.00 \times 10^1$ pg/ μ L

Limit of quantitation: 2.00×10^{-3} pg/ μ L

Precision: CV% \leq 15%

Standard curve



Datasheet

| Concentration (fg/ μ l) | Log ₁₀ Concentration | Ct Value(1) | Ct Value(2) | Ct Value(3) | Ct Mean Value | Recovery rate |
|-----------------------------|---------------------------------|-------------|-------------|-------------|---------------|---------------|
| 2.00E+04 | 4.30 | 15.30 | 15.23 | 15.21 | 15.24 | 0.30% |
| 2.00E+03 | 3.30 | 18.39 | 18.28 | 18.18 | 18.28 | 0.56% |
| 2.00E+02 | 2.30 | 21.86 | 21.82 | 21.78 | 21.82 | 0.18% |
| 2.00E+01 | 1.30 | 25.22 | 25.22 | 25.09 | 25.18 | 0.31% |
| 2.00E+00 | 0.30 | 28.55 | 28.58 | 28.43 | 28.52 | 0.29% |
| Amplification efficiency | | | | | | 99.1% |