

CHO Residual DNA Detection Kit (qPCR)

Overview

Cat.No. HG-CH001

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

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

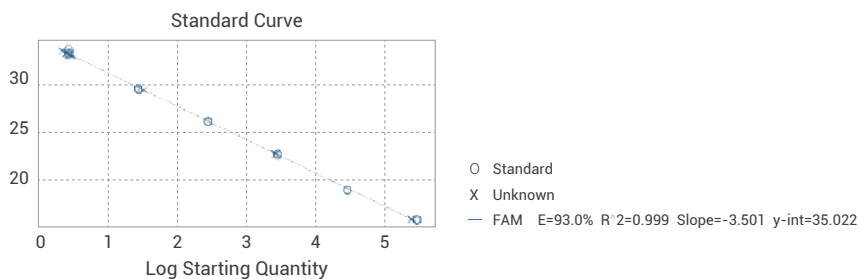
Specification

Assay range: 3.00~3.00x10⁵fg/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	15.94	16.08	16.12	16.05	0.59%
3.00E+04	4.48	19.10	19.16	19.17	19.14	0.18%
3.00E+03	3.48	22.72	22.89	22.95	22.86	0.51%
3.00E+02	2.48	26.26	26.38	26.30	26.31	0.23%
3.00E+01	1.48	29.61	29.70	29.79	29.70	0.30%
3.00E+00	0.48	33.88	33.28	33.43	33.53	0.93%
Amplification efficiency						93.0%