

Mycoplasma DNA Detection Kit (qPCR)

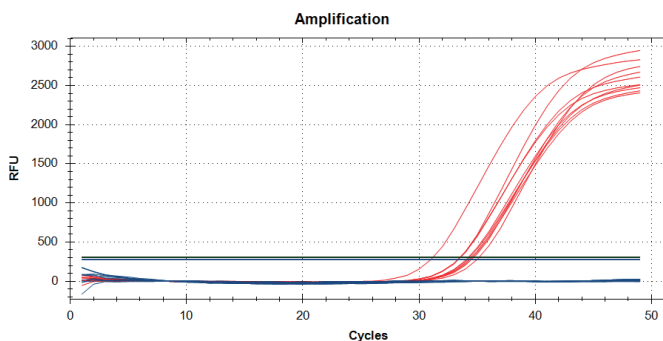
Overview

Cat.No. HG-ZY001

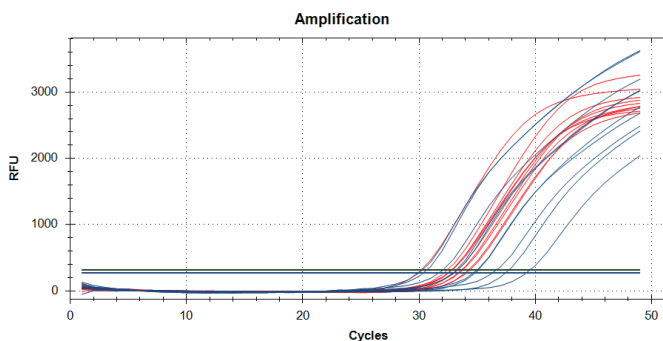
This kit is designed for the detection of mycoplasma contamination in master cell bank, working cell bank, cells for clinical use and biological products. This kit conforms to relevant regulations about mycoplasma testing in EP2.6.7 and JP XVI.

This kit adopts the qPCR-fluorescent probe method. The kit is a rapid, specific and reliable device and can finish the detection within 2 hours.

Test curve



Negative



Positive

— FAM
— HEX

Datasheet

	FAM(CT Value)		HEX(CT Value)	
	Negative	Positive	Negative	Positive
1	NA	31.72	34.45	32.62
2	NA	33.05	31.08	33.12
3	NA	36.32	33.43	33.45
4	NA	30.01	34.56	30.02
5	NA	30.37	35.05	33.55
6	NA	34.68	34.22	34.19
7	NA	34.83	34.35	32.63
8	NA	39.23	34.12	32.58
9	NA	32.95	33.39	34.07
10	NA	37.44	33.41	32.24

Results of 10 mycoplasma standards				Results of 3 relevant bacteria		
Strain	Positive/Total	Strain	Positive/Total	L.acidophilus	S.pneumoniae	Clos. acetobutyleum
M. orale	24/24	M. synoviae	24/24	Negative	Negative	Negative
M. gallisepticum	23/24	M. arginini	23/24			
A. laidlawii	24/24	M. hyorhinis	24/24			
M. fermentans	23/24	Spiroplasma citri	24/24			
M. pneumonia	24/24	M. saliarium	24/24			