

TaqMan® Pathogen Detection Kit

Reaction Setup

STEP	ACTION																																																																
1	<p>A. Make the Pathogen PCR Cocktail (●) as shown in this table.</p> <table border="1" style="margin: 10px 0;"> <thead> <tr> <th>Number of Unknowns</th> <th>TaqMan Pathogen Mix (μL)</th> <th>AmpliTaq Gold® DNA Polymerase (μL)</th> <th>Total (μL)</th> </tr> </thead> <tbody> <tr><td>10</td><td>880</td><td>20</td><td>900</td></tr> <tr><td>15</td><td>1144</td><td>26</td><td>1170</td></tr> <tr><td>20</td><td>1364</td><td>31</td><td>1395</td></tr> <tr><td>25</td><td>1628</td><td>37</td><td>1665</td></tr> <tr><td>30</td><td>1848</td><td>42</td><td>1890</td></tr> <tr><td>35</td><td>2112</td><td>48</td><td>2160</td></tr> <tr><td>40</td><td>2332</td><td>53</td><td>2385</td></tr> <tr><td>45</td><td>2596</td><td>59</td><td>2655</td></tr> <tr><td>50</td><td>2816</td><td>64</td><td>2880</td></tr> <tr><td>55</td><td>3080</td><td>70</td><td>3150</td></tr> <tr><td>60</td><td>3300</td><td>75</td><td>3375</td></tr> <tr><td>65</td><td>3564</td><td>81</td><td>3645</td></tr> <tr><td>70</td><td>3784</td><td>86</td><td>3870</td></tr> <tr><td>75</td><td>4048</td><td>92</td><td>4140</td></tr> <tr><td>80</td><td>4268</td><td>97</td><td>4365</td></tr> </tbody> </table> <p>B. Dispense 45 μL of the Pathogen PCR Cocktail into all of the wells except the NAC wells.</p> <p>This example shows 30 UNKN wells.</p>	Number of Unknowns	TaqMan Pathogen Mix (μL)	AmpliTaq Gold® DNA Polymerase (μL)	Total (μL)	10	880	20	900	15	1144	26	1170	20	1364	31	1395	25	1628	37	1665	30	1848	42	1890	35	2112	48	2160	40	2332	53	2385	45	2596	59	2655	50	2816	64	2880	55	3080	70	3150	60	3300	75	3375	65	3564	81	3645	70	3784	86	3870	75	4048	92	4140	80	4268	97	4365
Number of Unknowns	TaqMan Pathogen Mix (μL)	AmpliTaq Gold® DNA Polymerase (μL)	Total (μL)																																																														
10	880	20	900																																																														
15	1144	26	1170																																																														
20	1364	31	1395																																																														
25	1628	37	1665																																																														
30	1848	42	1890																																																														
35	2112	48	2160																																																														
40	2332	53	2385																																																														
45	2596	59	2655																																																														
50	2816	64	2880																																																														
55	3080	70	3150																																																														
60	3300	75	3375																																																														
65	3564	81	3645																																																														
70	3784	86	3870																																																														
75	4048	92	4140																																																														
80	4268	97	4365																																																														
2	<ul style="list-style-type: none"> • Add 5 μL Negative Control into each of the NTC wells. • Cap the 1st column of the plate with Optical caps. 																																																																
3	<ul style="list-style-type: none"> • Add 5 μL of your sample DNA into each of the UNKN (●) wells. • Cap all of the wells except those in column 12. 																																																																
4	<ul style="list-style-type: none"> • Add 44 μL TaqMan Pathogen Mix into the six NAC wells. 																																																																
5	<ul style="list-style-type: none"> • Add 5 μL Negative Control into the NAC wells. • Add 5 μL Positive Control into each of the two positive control wells. • Cap the wells in column 12. 																																																																

TaqMan® Pathogen Detection Kit

Procedure Flowchart

