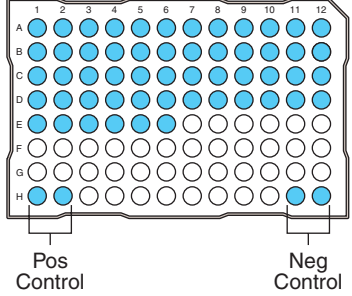
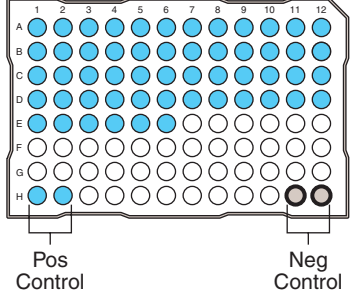
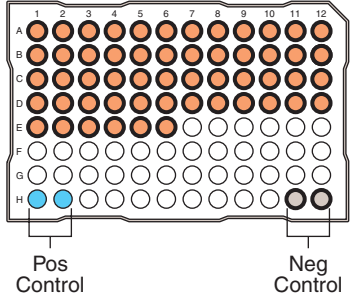
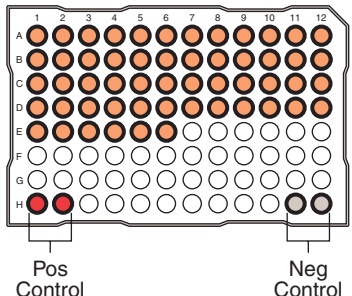


TaqMan® Pathogen Detection Kit

Reaction Setup

STEP	ACTION																																																																								
<h1>1</h1>	<p>⚠ WARNING BIOHAZARD. Biological samples such as tissues and blood have the potential to transmit infectious diseases. Read and follow the guidelines published in <i>Biosafety in Microbiological and Biomedical Laboratories</i> (http://bmlb.od.nih.gov) and in the Occupational Safety and Health Standards, Bloodborne pathogens (http://www.access.gpo.gov/nara/cfr/waisidx_01/29cfr1910a_01.html). Wear appropriate protective eyewear, clothing, and gloves.</p> <table border="1" data-bbox="267 436 722 850"> <thead> <tr> <th>Number of Reactions</th> <th>Pathogen PCR Mix (μL)</th> <th>AmpliTaq Gold® DNA Polymerase (μL)</th> <th>Total Volume</th> </tr> </thead> <tbody> <tr><td>10</td><td>484</td><td>11</td><td>495</td></tr> <tr><td>15</td><td>748</td><td>17</td><td>765</td></tr> <tr><td>20</td><td>968</td><td>22</td><td>990</td></tr> <tr><td>25</td><td>1188</td><td>27</td><td>1215</td></tr> <tr><td>30</td><td>1452</td><td>33</td><td>1485</td></tr> <tr><td>35</td><td>1672</td><td>38</td><td>1710</td></tr> <tr><td>40</td><td>1936</td><td>44</td><td>1980</td></tr> <tr><td>45</td><td>2156</td><td>49</td><td>2205</td></tr> <tr><td>50</td><td>2376</td><td>54</td><td>2430</td></tr> <tr><td>55</td><td>2640</td><td>60</td><td>2700</td></tr> <tr><td>60</td><td>2860</td><td>65</td><td>2925</td></tr> <tr><td>65</td><td>3124</td><td>71</td><td>3195</td></tr> <tr><td>70</td><td>3344</td><td>76</td><td>3420</td></tr> <tr><td>75</td><td>3564</td><td>81</td><td>3645</td></tr> <tr><td>80</td><td>3828</td><td>87</td><td>3915</td></tr> <tr><td>85</td><td>4048</td><td>92</td><td>4140</td></tr> <tr><td>90</td><td>4312</td><td>98</td><td>4410</td></tr> </tbody> </table> <p>A. Make the pathogen PCR Cocktail (●) as shown in this table.</p> <p>B. Dispense 45 μL of the pathogen PCR Cocktail into all of the designated wells.</p> <p>This example shows the plate setup for qualitative (pathogen presence or absence) analysis mode.</p> <p>For quantitative analysis mode, refer to the User Guide, Chapters 3, 4, and 6.</p> 	Number of Reactions	Pathogen PCR Mix (μL)	AmpliTaq Gold® DNA Polymerase (μL)	Total Volume	10	484	11	495	15	748	17	765	20	968	22	990	25	1188	27	1215	30	1452	33	1485	35	1672	38	1710	40	1936	44	1980	45	2156	49	2205	50	2376	54	2430	55	2640	60	2700	60	2860	65	2925	65	3124	71	3195	70	3344	76	3420	75	3564	81	3645	80	3828	87	3915	85	4048	92	4140	90	4312	98	4410
Number of Reactions	Pathogen PCR Mix (μL)	AmpliTaq Gold® DNA Polymerase (μL)	Total Volume																																																																						
10	484	11	495																																																																						
15	748	17	765																																																																						
20	968	22	990																																																																						
25	1188	27	1215																																																																						
30	1452	33	1485																																																																						
35	1672	38	1710																																																																						
40	1936	44	1980																																																																						
45	2156	49	2205																																																																						
50	2376	54	2430																																																																						
55	2640	60	2700																																																																						
60	2860	65	2925																																																																						
65	3124	71	3195																																																																						
70	3344	76	3420																																																																						
75	3564	81	3645																																																																						
80	3828	87	3915																																																																						
85	4048	92	4140																																																																						
90	4312	98	4410																																																																						
<h1>2</h1>	<ul style="list-style-type: none"> • Add 5 μL Negative Control into each of the two negative control wells. • Cap the wells. 																																																																								
<h1>3</h1>	<ul style="list-style-type: none"> • Add 5 μL sample DNA into each of the designated UNKN (●) wells. Note: This example shows 54 sample reactions with four control reactions. • Cap the wells. 																																																																								
<h1>4</h1>	<ul style="list-style-type: none"> • Add 5 μL Positive Control into each of the two positive control wells. • Cap the wells. 																																																																								

TaqMan® Pathogen Detection Kit

Procedure Flowchart

