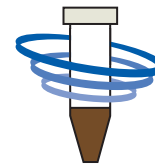
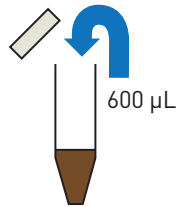
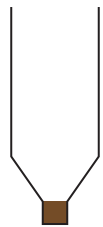


Pathatrix® system—VIDAS® strip linking step

Standard sample



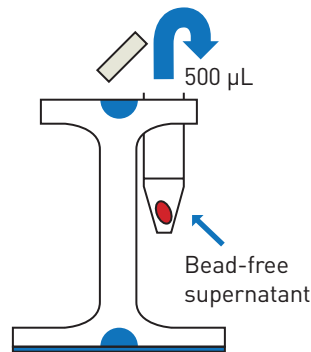
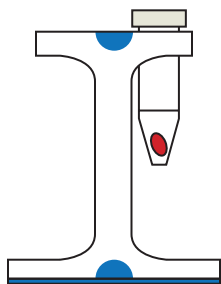
Add **700 µL** of **pre-warmed BPW (37°C)** to resuspend the Pathatrix® beads

Transfer **600 µL** of bead suspension into a standard 2.0 mL screw cap microfuge tube

Vortex the tube for 10 sec and then incubate the bead suspension for **2 hr at 37°C**

Vortex the tube for 10 sec and then place in a water bath at **95°C** and heat for **15 min** to release antigens

(Note: 100 µL of the beads can be retained for plating)



Immediately after heating, capture beads for **1 min** on Magnetic Capture Rack

Remove **500 µL** of **bead-free supernatant** and inoculate the VIDAS® strip

Proceed with the assay in accordance with the manufacturer's instructions