

# Pathatrix® system workflow—sample preparation and loading



1. Weigh samples into pre-warmed pre-enrichment broth.

Incubate the samples at the correct temperature and time.



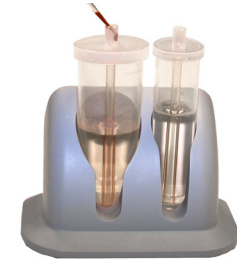
2. Prepare sample pools using Pooling Straws and Filta-Foam.

Place your sample (**65 mL maximum**) into the Sample Vessel.



3. Pour sterile PBS into the Elution Vessel to the fill line (35 mL).

Push the appropriate lids firmly onto the vessels.



4. Add a 50 µL bead dose into the Sample Vessel lid spout.

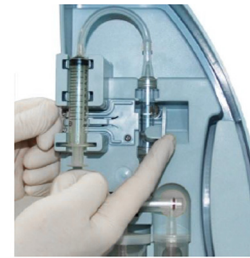


5. Remove the sterile kit from the bag and orient so the plunger points left.

Place the valve into the lid spouts and push down firmly.



6. Carefully holding the vessels to avoid spills, lift the vessels and attached consumable and insert the vessels into the cartridge, ensuring the base of the Elution Vessel is fully pushed back. Working upward, push the phase into the cartridge and put the syringe in place.



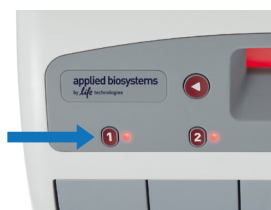
7. Slide the magnet assembly into locking position and check that it will release using the button on the side of the cartridge.

Re-slide the magnet assembly across into the locking position.



8. Insert the cartridge into the Pathatrix® system until locked into place.

# Pathatrix<sup>®</sup> system workflow—running test and bead isolation



1. Press the numbered button above the loaded cartridge to begin the run.



2. Once the run step has finished, the LED will flash **green** and **red**. Press the numbered button to initiate the draining step.



3. When the draining step is complete, the LED will turn **red**. Pull the cartridge firmly out of the instrument.



4. Place the cartridge in the rack provided and remove the consumable and vessels, starting with the syringe and working downward.

Hold onto the Sample and Elution Vessels to prevent spillage.



5. Place the vessels into the Tube Rack.

Detach the lid from the Elution Vessel and discard the consumable and attached Sample Vessel.



6. Leave the Elution Vessel in the rack for 1 min to allow the beads to be collected on the magnet.



7. Without removing the Elution Vessel from the rack, remove all of the wash buffer from the Elution Vessel, leaving the beads in place.



8. Resuspend the beads using 100  $\mu$ L of PBS, and proceed to detection methodology.