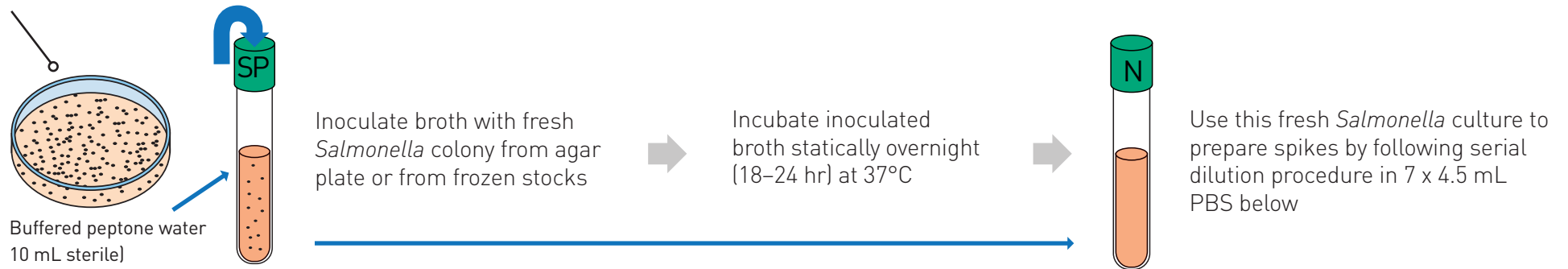
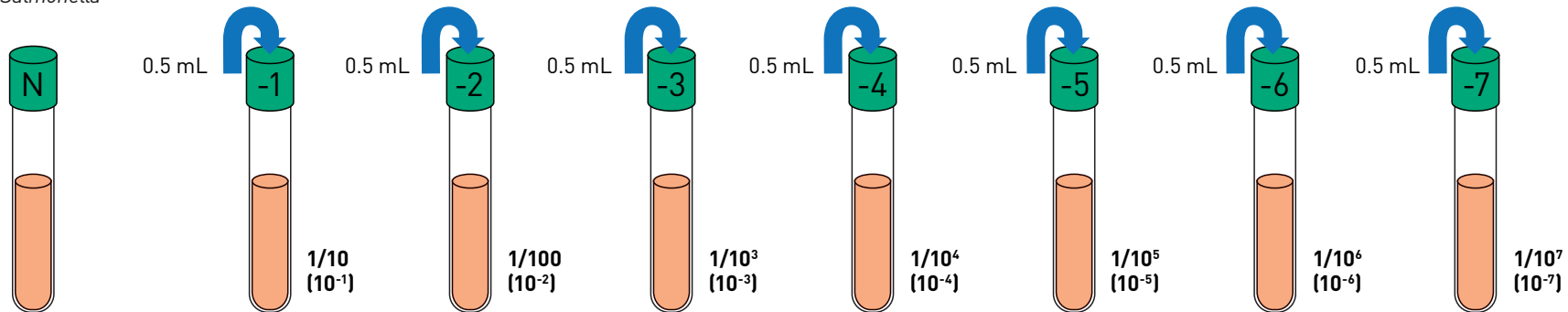


Creating a *Salmonella* culture serial dilution



Overnight culture of *Salmonella*

7 sterile tubes (e.g., plastic bijoux) each containing 4.5 mL sterile PBS



Starting with the fresh overnight *Salmonella* culture and using sterile technique, pipette 0.5 mL from each tube into the next sterile PBS tube. After liquid transfer, secure cap and mix well before taking next 0.5 mL. Change pipette tip between transfers.

Plate 2 x 100 µL from dilutions 6 and 7 only onto 2 well-dried agar plates (e.g., XLD or TSA)

Incubate plates and count colonies next day

