

# Attention

**Please do not remove the drainage tubing once installed. Removal of the drainage tubing may cause the drain port fittings to loosen, leading to leaks. Cleaning can be easily performed with the tubing attached. Please see the instruction manual for additional maintenance procedures.**

## **Assembling the drainage system:**

1. Remove the stopcocks and tubing from the bag.
2. Slide the white clamps over the short pieces of tubing (with the stopcocks already attached). Don't tighten them yet.
3. Attach the ends of the short pieces of tubing to both drain port fittings at the base of the solution tank.

**Important:** Do not attach the tubing by twisting or rotating in a circular motion, this may loosen the drain port fittings and cause leaks when the tank is filled with solution. Instead bend the tubing up and down while applying pressure to push the tubing onto the fittings.

4. Now, tighten the white clamps to ensure a complete seal and to prevent leaks.

**Note:** It may be easier to use pliers to tighten the white clamps.

5. Attach the long pieces of drainage tubing to the other end of the stopcocks. Ensure the tubing reaches your desired receptacle/ reservoir.
6. Make sure the stopcocks are in the closed position prior to pouring in any solutions.