

Instructions for Assembling OmicsArray™ Slide Incubation Chamber Set (PA105)

1. Place the glass slide over the base, ensure the array side is facing up (the line mark should be on the bottom left side as show in the figure 1).



Figure 1

2. Place the chamber and gasket module on the slide, as shown in the figure 2.



Figure 2

Attention: When placing the chamber and gasket module, it must be placed in the groove and kept flat, otherwise it may cause the damage of the array.

3. Snap the side clips as shown in Figure 3.

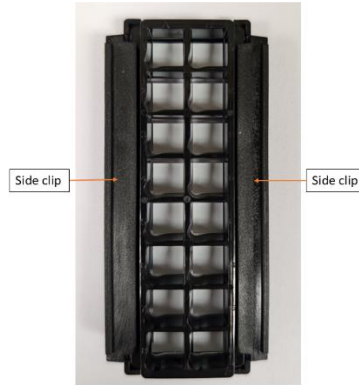


Figure 3

4. Install all 3 chamber modules according to the instruction step 1-3 and place the assembled modules in chamber rack (Figure 4).



Figure 4

5. Cover the chamber with the lid, as shown in figure 5.



Figure 5