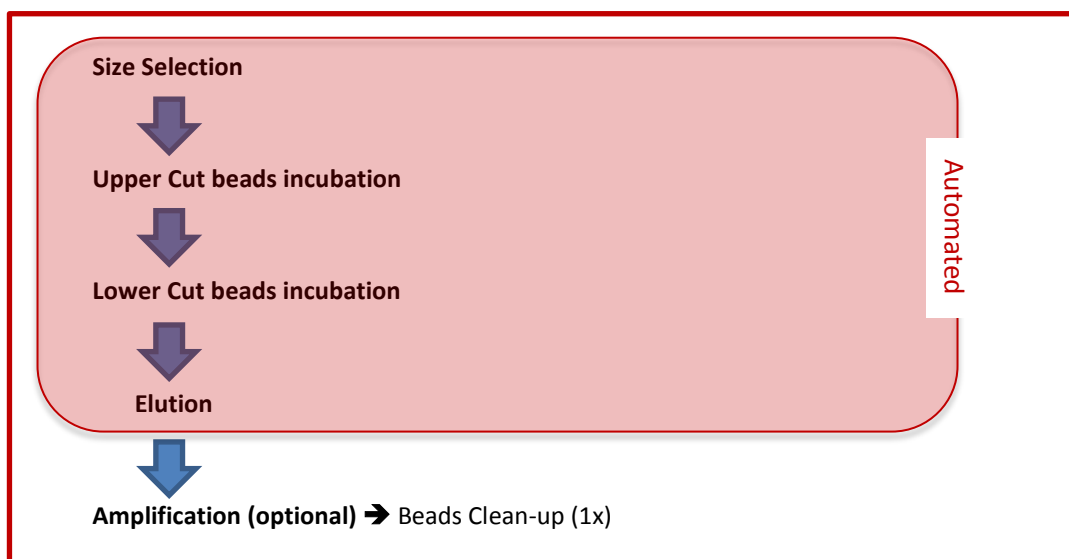


1. About the protocol

The “AMPure_XP_Size_Selection” protocol on the IP-Star[®] is using the standard “Agencourt[®] AMPure[®] XP Beads” from Agencourt[®]. It allows you to perform size selection from 250bp to 500bp just with changing the amount of beads.

It provides flexibility to perform 1 to 16 samples in one run starting with **100 µl** of sample. The whole protocol takes approximately 1h for 8 samples. At the end, you recover size selected products ready for amplification.

2. Workflow



3. Material required

a. Reagents & kits

Item	Supplier
Agencourt [®] AMPure [®] XP Beads	Beckman Coulter [®]
Fresh Ethanol 80%	Lab supplier

b. Consumables

Item	Supplier	Catalogue #
200 µl tube strips (8 tubes/strip) + cap strips	Diagenode	C30020002
2 ml microtube	Diagenode	C30010014
Medium reagent container	Diagenode	C30020003
96 well microplates	Diagenode	C30080030
Tips (box)	Diagenode	C30040021
Tips (bulk)	Diagenode	C30040020

4. IP-Star setup

- Switch ON the IP-Star.
- Select “**Protocols**” icon and then click on “**Library prep**”.
- Under “**Library prep**”, select “**AMPure_XP_Size_Selection**”.

Note:

If you plan to run between 1 and 8 samples, chose “**AMPure_XP_Size_Selection_08**”

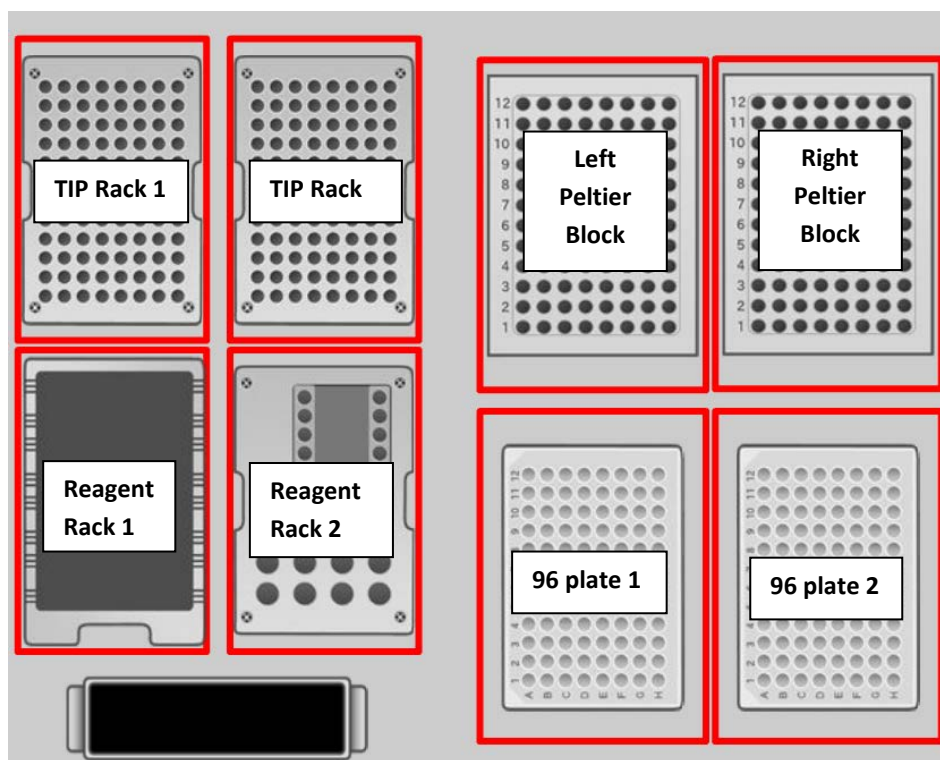
If you plan to run between 9 and 16 samples, chose “**AMPure_XP_Size_Selection_16**”

- Setup the exact number of samples that you want to process.

Note:

The **Left Peltier Block** is now cooling down to 4°C to keep your samples cold.

- Setup all the plastics on the platform according to the screen layout.



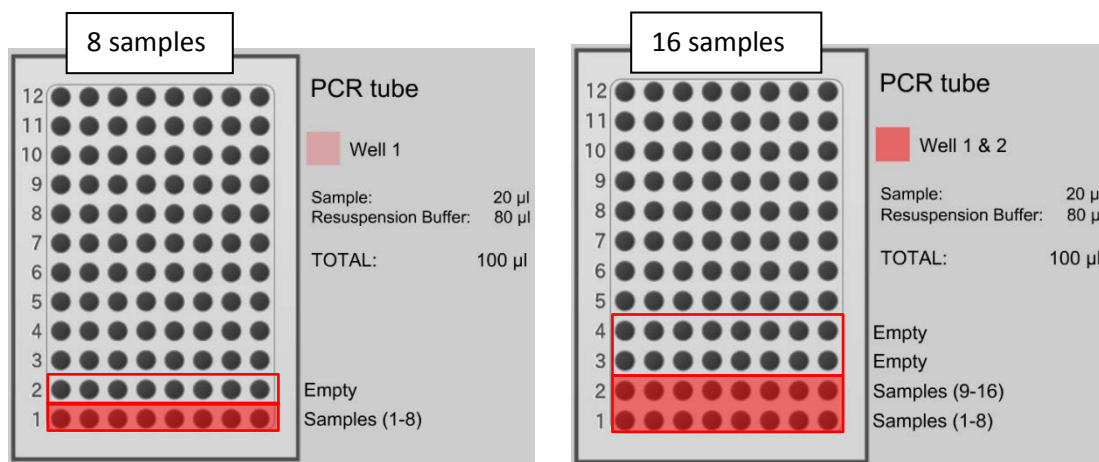
- Fill **TIP Rack 1** (and 2 if processing 16 protocol) with tips according to the screen.
- Fill **Reagent Rack 1 & 2** with reagent containers according to the screen.
- Fill **96 plate 1** with 96 well microplates.
- Fill **Left Peltier Blocks** with 200 µl tube strips according to the screen.

5. Reagents & Samples setup

Note:

Allow “Agencourt[®] AMPure[®] XP Beads” to come at room temperature.

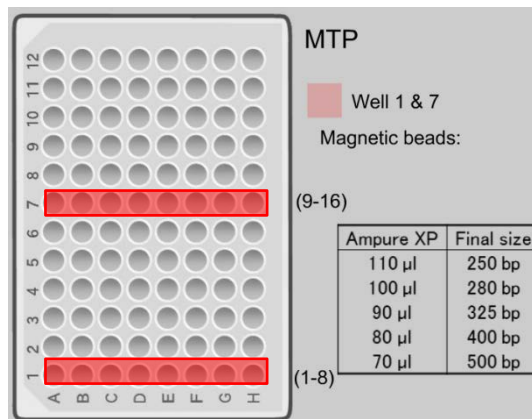
- Fill the **Left Peltier Block** with the samples according to the screen layout.
- Fill **20 µl of Samples** in lane 1 (and 2 if processing more than 8 samples).
- Fill up to 100µl with 80µl of resuspension buffer.



- Fill **Agencourt[®] AMPure[®] XP Beads** in lane 1 on **96 Plate 1** (and 7 if processing more than 8 samples) according to the required size following recommendations from the table.

Note:

Resuspend the beads with pipetting up and down several times before dispense them.



- Fill **freshly prepared Ethanol 80%** in the container on the **Reagent Rack 1**.
- Fill **Resuspension Buffer** in the container on **Reagent Rack 2**.
- Close the door and Run.

6. End

- Recover your samples on the **Left Peltier Block**