



Item No.: H1010-MR

Thank you for purchasing the Incu-Shaker 10L with MAGic Clamp Platform.

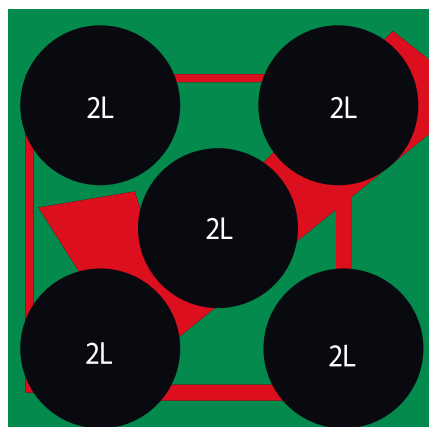
For your convenience, the MAGic Clamp platform has already been installed to your incubator.

Flask Clamp Installation and Alignment

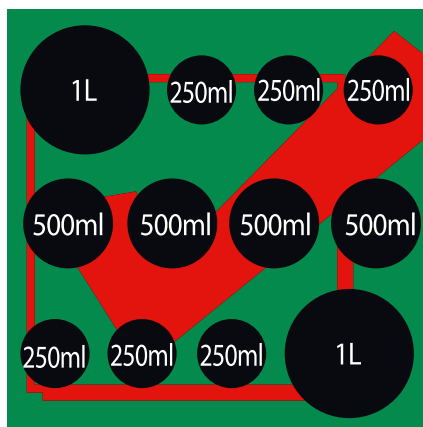
The MAGic Clamp Platform is capable of supporting 5 x 2000ml, 9 x 1000ml, 14 x 500ml, 25 x 250ml 36 x 125ml, 49 x 50ml flasks or various combinations of these flask sizes.

In addition, this platform offers several accessories racks for securing tubes and bottle.

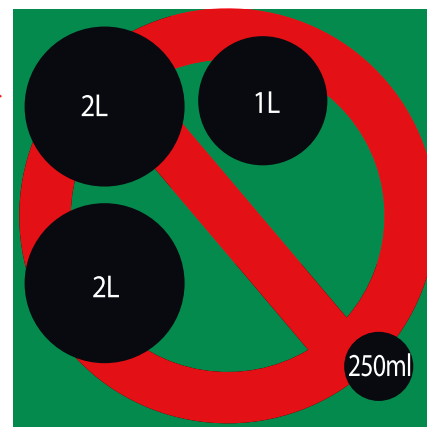
To attach a flask or tube rack clamp, simply place the clamp (magnet side down) on the platform in the desired location. When attaching clamps, **always** ensure a balanced load. (Please see below for an example of both properly and improperly balanced loads.



Example of a properly balanced load containing 5x2L flasks



Example of properly a balanced load containing a mixture of flask sizes



Example of an improperly a balanced load

Benchmark 
Scientific

TEL: 908-769-5555 FAX: 908-222-1864 EM: info@BenchmarkScientific.com WEB: www.BenchmarkScientific.com