

## Capit-All and Capit-All IS Screw Cap Tube Cappers/Decappers

Thermo Scientific and The Automation Partnership (TAP) Capit-All and Capit-All IS instruments remove and replace screw caps on 96- and 24-format storage tubes with precision and ease. The instruments' robust designs ensure reliable, high throughput access and storage of samples in Thermo Scientific Matrix, NUNC and other screw cap storage tubes, while reducing risk of repetitive strain injury.



**Enhanced Productivity.** Increase laboratory throughput and improve ergonomics by capping or decapping up to 96 tubes at once (in approximately 10 seconds/rack). With the addition of spacers, enable use with 500  $\mu$ l or 1.0 ml tube racks on the same instrument and maintain seal integrity with pre-determined torque control.

**Superior Sample Protection.** Safety features ensure that racks are loaded in the correct orientation, as well as prevent cross-threading or overtightening of caps. A drip tray prevents cross contamination of samples during entry and removal. The instruments are also engineered to be cleaned and maintained without disassembly to further ensure samples remain debris free.

**Ease of Use.** The simple, four-button user interface enables laboratory personnel of any level to operate the instruments with minimal training required.

**Ideal for Scale up.** The Capit-All<sup>®</sup> IS can be operated as a stand-alone benchtop unit. When research needs to be ramped up, the instrument easily can be integrated with most any robotic system. The Capit-All IS can also process over one million tubes before routine service to maximize uptime.

**Knowledgeable Support.** Our team has a strong reputation for supporting both equipment and consumables, and can demonstrate simple ways to improve sample management and productivity in your laboratory using the Capit-All or Capit-All IS Screw Cap Tube Capper/Decapper.

## Capit-All and Capit-All IS Screw Cap Tube Cappers/Decappers

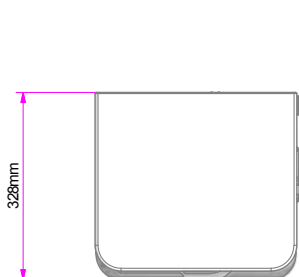
	Capit-All	Capit-All IS
Dimensions (W x D x H)	12.9"/360 mm x 14.2"/328 mm x 22.0"/560 mm	27.6"/700 mm x 13.4"/340 mm x 22.0"/560 mm
Weight	97 lbs/45 kg	129 lbs/58.5 kg
Power Supply	110-240 V AC, 50-60 Hz , 6 Amps	110-240 V AC, 50-60 Hz, 6 Amps
Speed	Approx. 10 seconds per cycle (cap or decap)	Approx. 15 seconds per cycle (cap or decap)
Operation	Hand operation via push buttons	Remotely via RS232; By computer via included drivers; Hand operation via push buttons

### Ordering Guide

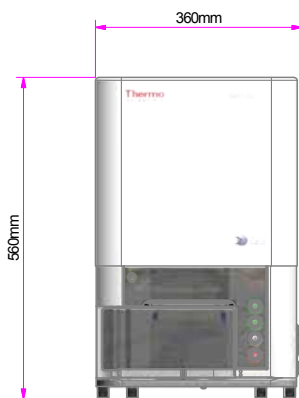
Item No.	Description
4111MAT	Capit-All 96-Channel Screw Cap Tube Capper/Decapper for use with Matrix® style screw top caps
4111NUN	Capit-All 96-Channel Screw Cap Tube Capper/Decapper for use with Nunc® style screw top caps
4112NUN	Capit-All 24-Channel Screw Cap Tube Capper/Decapper for use with Nunc style internal thread cryostorage tubes
4113NUN	Capit-All 24-Channel Screw Cap Tube Capper/Decapper for use with Nunc style external thread cryostorage tubes
4113NAL	Capit-All 24-Channel Screw Cap Tube Capper/Decapper for use with Nalgene® style external thread cryostorage tubes
4111MAT-IS	Capit-All IS 96-Channel Automation Friendly Screw Cap Tube Capper/Decapper for use with Matrix style screw top caps
4111NUN-IS	Capit-All IS 96-Channel Automation Friendly Screw Cap Tube Capper/Decapper for use with Nunc style screw top caps

If a model is required for a different style of cap, please speak with your sales representative

#### Capit-All:

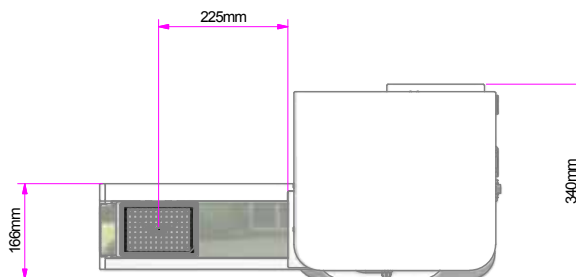


▲ Figure 1: Top View

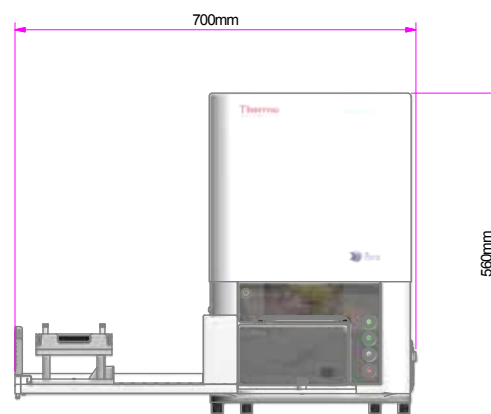


▲ Figure 1: Front View

#### Capit-All IS:



▲ Figure 2: Top View



▲ Figure 2: Front View