

## Product Guidelines for Site Installation

Product Name:	Thermo Scientific™ Barnstead™ Mega-Pure® Water Distillation Systems	Item Description:	Tap to Type 2 Lab Water System
Model #	MP-3A, 3.4LPH	Manufacturer	Thermo Scientific
Order #	A440367, A440696		

### A. RECOMMENDED FEED WATER SPECIFICATIONS

Resistivity	µS/cm	> 0.001	µS/cm
Blocking Index	SDI max.	<3	SDI
Manganese	mg/l	> 0.05	mg/l
Iron	mg/l	> 0.05	mg/l
Free Chlorine	mg/l	< 0.1	mg/l
pH Range	pH	4-11	pH
Pressure	PSI	20-100	PSI
Flow Rate	GPH	4	GPH
Recommended Type	ASTM 1, 2, or 3	Tap	DIN2000

### B. ELECTRICAL POWER REQUIREMENTS

Electrical Specifications								
<b>A440696</b>	208	Volts	50/60	Hertz	13	Amps	Single	# of Phases
<b>A440367</b>	240	Volts	50/60	Hertz	11	Amps	Single	# of Phases
Total number of receptacles / plugs required to operate water system.							1	
Length of line cord	5	Feet	0	Inches				

### C. PLUMBING REQUIREMENTS

Water Inlet Line	
Type of Fitting	Male NPT
Size of Fitting	1/8
Manual Cutoff Valve	Strongly Recommend
Waste Drain	
Type of Drain	Atmospheric
Max. Distance from Unit	5
	Feet

### D. SPATIAL REQUIREMENTS

Equipment exterior dimensions (inches):	23	Width	12	Depth	45	Height
Automatic Collection System (ACS) Dimensions (inches):	16.25	Width	16.25	Depth	38	Height
Dual Cartridge Deionizer (D2) Dimensions (inches):	16.75	Width	7.5	Depth	25.5	Height
Equipment Weight (lbs):	60	Lbs.				
ACS Reservoir Weight (full) (lbs):	60	Lbs.				
Dual Cartridge Deionizer Weight (lbs):	30	Lbs.				

Wall mount specifications: If the **system** is to be mounted on the wall or on a shelf the designated area must be capable of supporting up to 200 lbs.

### E. Accessories / Replacement Consumables

#### Required Accessories

#### Reservoir Select One from the three options below:

- 413964 6L Plastic Reservoir
- 410535 9L Glass Reservoir
- 413934 13L Glass Reservoir
- 410164 45L Glass Reservoir
- D440704 Automatic Collection System (includes 45L glass reservoir)
- 4402362 Dual Solenoid required if using pretreated boiler feed water and untreated cooling water

#### Optional Accessories

- D440066 Dual Cartridge Deionizer (**REQUIRES** Still Adapter Kit (#440376))

#### Cartridges Select 2 from the three options below:

- D400377 High Purity Cartridge
- D400499 High Capacity Cartridge
- D440265 Organic Removal Cartridge