

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-6A, 6LPH	Manufacturer	Thermo Scientific
Order #	A440518, A440697		

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								
A440696	208	Volts	50/60	Hertz	25	Amps	Single	# of Phases
A440367	240	Volts	50/60	Hertz	21	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 Inch
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:

- 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