

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-11A, 13LPH	Manufacturer	Thermo Scientific
Order #	A440118, A440117		

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								
A440117	208	Volts	50/60	Hertz	49	Amps	Single	# of Phases
A440118	240	Volts	50/60	Hertz	42	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):	29.2	Width	14.4	Depth	53	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