

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-12A, 12LPH	Manufacturer	Thermo Scientific
Order #	A442011, A442012		

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							
A442012	208	Volts	50/60	Hertz	49	Amps	Single # of Phases
A442011	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)

Optional Accessories

400634	Automatic Collection System Wall Mounting Bracket
--------	---