

Product Guidelines for Site Installation

Product Name:	Thermo Scientific™ Barnstead™ Classic Water Distillation Systems	Item Description:	Tap to Type 2 Lab Water System
Model #	5 GPH	Manufacturer	Thermo Scientific
Order #	A1015-C, A1015-B-61		

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	40-90	PSI	
Flow Rate	GPH	4	GPH	
Recommended Type	ASTM 1, 2, or 3	Tap	DIN2000	

B. ELECTRICAL POWER REQUIREMENTS

Electrical Specifications								
A1015-C	208	Volts	50/60	Hertz	36	Amps	Single	# of Phases
A1015-B-61	240	Volts	50/60	Hertz	57	Amps	Single	# of Phases
Total number of receptacles / plugs required to operate water system.							2	
Length of line cord		5	Feet	0	Inches			

C. PLUMBING REQUIREMENTS

Water Inlet Line	
Type of Fitting	Male NPT
Size of Fitting	1/4 Inch
Manual Cutoff Valve	Strongly Recommend
Waste Drain	
Type of Drain	Atmospheric
Max. Distance from Unit	3 Feet
Size of Drain Fitting	1/2"
The drain construction should be able to handle 180°C water and 40 GPH	

D. SPATIAL REQUIREMENTS

Equipment exterior dimensions (inches):	35	Width	14	Depth	77	Height
Equipment Weight (lbs):	80	Lbs.				

E. Accessories / Replacement Consumables

Required Accessories

H3111	Ventgard Air Filter and Water Seal
H4005	UV Light Assembly

Reservoir **Select One from the options below:**

B3043	10 Gallon Tin Lined Reservoir
B3045	25 Gallon Tin Lined Reservoir
B3046	50 Gallon Tin Lined Reservoir
B3047	100 Gallon Tin Lined Reservoir
B3049	200 Gallon Tin Lined Reservoir

Reservoir Floor Stand* or Wall Mounting Bracket (Select One from the options below):

H1000	Floor Stand - 10 Gallon Reservoir
H1001	Floor Stand - 25 Gallon Reservoir
H1002	Floor Stand - 50 Gallon Reservoir
H1003	Floor Stand - 100 Gallon Reservoir
H3230	Floor Stand - 200 Gallon Reservoir
H3240	Wall Mounting Bracket - 10 Gallon Reservoir
H3242	Wall Mounting Bracket - 25 Gallon Reservoir

* Customers are responsible for securing stand to floor.

Optional Accessories

G2110	Fully Automatic Controls
-------	--------------------------