

## Product Guidelines for Site Installation

Product Name: **Thermo Scientific™ Barnstead™ Classic Water Distillation Systems** Item Description: **Tap to Type 2 Lab Water System**

Model # **1 GPH** Manufacturer **Thermo Scientific**  
Order # **A1011-A, A1011-B-61**

### A. RECOMMENDED FEED WATER SPECIFICATIONS

Resistivity	µS/cm	<b>&gt; 0.001</b>	µS/cm
Blocking Index	SDI max.	<b>&lt;3</b>	SDI
Manganese	mg/l	<b>&gt; 0.05</b>	mg/l
Iron	mg/l	<b>&gt; 0.05</b>	mg/l
Free Chlorine	mg/l	<b>&lt; 0.1</b>	mg/l
pH Range	pH	<b>4-11</b>	pH
Pressure	PSI	<b>40-90</b>	PSI
Flow Rate	GPH	<b>4</b>	GPH
Recommended Type	ASTM 1, 2, or 3	<b>Tap</b>	DIN2000

### B. ELECTRICAL POWER REQUIREMENTS

Electrical Specifications							
<b>A1011-A</b>	<b>120</b>	Volts	<b>50/60</b>	Hertz	<b>23</b>	Amps	<b>Single</b> # of Phases
<b>A1011-B-61</b>	<b>240</b>	Volts	<b>50/60</b>	Hertz	<b>12</b>	Amps	<b>Single</b> # of Phases
Total number of receptacles / plugs required to operate water system.						<b>2</b>	
Length of line cord	<b>5</b>	Feet	<b>0</b>	Inches			

### C. PLUMBING REQUIREMENTS

Water Inlet Line			
Type of Fitting	<b>Male NPT</b>		
Size of Fitting	<b>1/4</b>	Inch	
Manual Cutoff Valve	<b>Strongly Recommend</b>		
Waste Drain			
Type of Drain	<b>Atmospheric</b>		
Max. Distance from Unit	<b>3</b>	Feet	
Size of Drain Fitting	<b>1/2"</b>		The drain construction should be able to handle 180°C water and 8 GPH

### D. SPATIAL REQUIREMENTS

Equipment exterior dimensions (inches):	<b>22</b>	Width	<b>10</b>	Depth	<b>66</b>	Height
Equipment Weight (lbs):	<b>60</b>	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

G2100	Fully Automatic Controls
-------	--------------------------