

## Product Guidelines for Site Installation

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

Model # **10 GPH** Manufacturer **Thermo Scientific**  
Order # **A1016-D**

### 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  
**A1016-D** **240** Volts **50/60** Hertz **66** 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 80 GPH

### D. SPATIAL REQUIREMENTS

Equipment exterior dimensions (inches): **37** Width **14** Depth **89** Height  
 Equipment Weight (lbs): **100** 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

G2125 Fully Automatic Controls