

TURBO-V HEATER BAND

Model 969-9801 to 969-9814

Model 969-9317, 969-9327



To reach base pressure in the chamber or system in the shortest possible time, it is recommended to bake out the chamber and at the same time to bake out the Turbo-V pump.

For this purpose an heater band is provided.

The heater is shaped to fit the upper part of the envelope (fig. 1) and will automatically heat it to a temperature of about 80°C.

The required bake out time will be a function of the degree of contamination of the pump and of the desired base pressure value.

Whenever possible it is recommended to connect the heater band using the facilities already built into the Turbo-V controller; otherwise plug it to the appropriate outlet socket.

 CAUTION

The maximum permissible temperature of the Turbo-V pump inlet flange is 120°C (250°F). If the system chamber is to be baked out at high temperature, a radiation shield must be used to prevent overheating of the inlet flange on the Turbo-V pump.

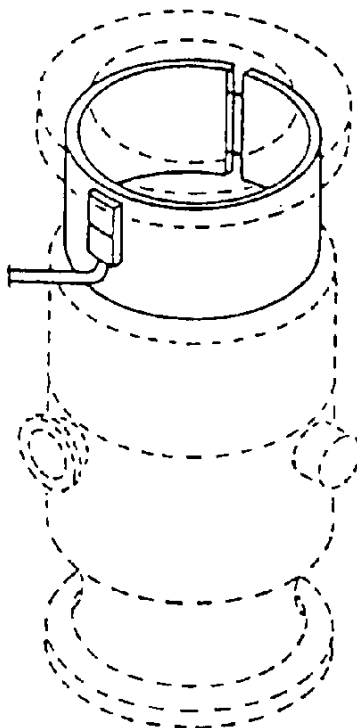


Fig. 1
Heater installation

The table reports the outline dimensions, specifications and the applicability of the heater bands.

HEATER BANDS	USED ON	INPUT VOLTAGE POWER		INTERNAL DIAMETER	HEIGHT
969-9801 969-9802	Turbo-V 70 series	220V 120V	50W 50W	80 (3.15)	20 (.787)
969-9801 969-9802	Turbo-V 80 series	220V 120V	50W 50W	80 (3.15)	20 (.787)
969-9803 969-9804	Turbo-V 200 series	220V 120V	60W 60W	116 (4.56)	30 (1.18)
969-9803 969-9804	Turbo-V 250 series	220V 120V	60W 60W	116 (4.56)	22 (.866)
969-9805 969-9806	Turbo-V 300 series	220V 120V	100W 100W	157 (6.18)	35 (1.37)
969-9803 969-9804	Turbo-V 300HT series	220V 120V	60W 60W	116 (4.56)	22 (.866)
969-9807 969-9808	Turbo-V 450 series	220V 120V	100W 100W	190 (7.48)	50 (1.96)
969-9807 969-9808	Turbo-V 500 series	220V 120V	100W 100W	190 (7.48)	50 (1.96)
969-9807 969-9808	Turbo-V 700HT series	220V 120V	100W 100W	190 (7.48)	50 (1.96)
969-9809 969-9310	Turbo-V 1000 series	220V 120V	200W 200W	234 (9.21)	70 (2.75)
969-9317 969-9327	Turbo-V 1000HT series	220V 120V	200W 200W	234 (9.21)	30 (1.18)
969-9811 969-9812	Turbo-V 1800 series	220V 120V	200W 200W	295 (11.61)	90 (3.54)
969-9813 969-9814	Turbo-V 3500 series	220V 120V	250W 250W	444 (17.48)	100 (3.93)

Dimensions mm
 inches