

SERVICE "S" PARAMETERS		
Accessing Service Mode: Set dial on remote to 0 and then hold the Receptacle button for 10 seconds		
Exiting Service Mode: Press and release Light button		
If unable to enter Service Mode check that the unit is not in OFF mode by pressing the Set button for 10 sec		
S 1	Downflow blower	0-100
S 2	Exhaust blower	0-100
S 3	Downflow alarm	0-100
S 5	Inflow alarm	0-100
S 7	Operating hours	Variable
S 8	Reset Operating Hours	Variable
S 9	UV hours / reset	Variable
S 10	Displayed DF	75
S 11	Displayed IF	105
S 12	Remote Address	0 - F
S 13	KI Discus	5
S 14	Time before alarm	2
S15	Window position code	See table below
S 16	Autocal	0-Off, 1-Short, 2-Long

Window Position Code		
Status front window	Indication Code	Check switch
Front window in maximal opening position	0100	Function is correct
	1100	S6 and/or S11
	0101	S7
Moving front window	1101	S7 and S6 and/or S11
	0000	Function is correct
	1000	S6, S11
Front window in working position	0001	S7
	1001	S7, S6, S11
	0010	Function is correct
	1010	S6, S11
	0011	S7
	0000	S2,S3
	0001	S2,S3,S7
Front window closed	1000	S2,S3,S6,S11
	1001	S2,S3,S7,S6,S11
	1011	S7, S6,S11
	0011	Function is correct
	0001	S2,S3
	0010	S7
	1011	S6, S11
	1010	S7, S6,S11
	1000	S2,S3,S7,S6,S11
	1001	S2,S3,S6,S11
	0000	S2,S3,S7

Operating Modes
For full description see page 6-3 of operators manual
Flourecent light will function in all modes. Outlet all but UV mode
OFF Mode - Blowers off. Display shows time.
Standby Mode - Blowers reduced. Display shows time.
Work Mode - Blowers at full speed.
UV Mode - Blowers reduced. Display shows countdown timer.

To enter the UV Timer press and hold the UV button until the ready signal sounds. The last run time that was selected will be displayed. Press up or down to adjust the time in 30min increments and then press Ctrl to save.

PROGRAM "P" PARAMETERS		
Accessing Parameter Mode: Hold Light button for 10 seconds while in Service mode		
Exiting Parameter Mode: Press and release Light button.		
P1	Software Version	2.1
P2	Parameter ID	See config table
P3	Window Type	0
P4	UV Lamp Socket (0-No 1-Yes)	0
P5	Potential Contacts	0-No / 1-Yes
P6	Monitor Alarm	0-No / 1-Yes
P7	Audible Alarm Silence	0-No/ 1-Yes/ 2-Dis
P8	Unit of Measure	0-m/s / 1-fpm
P9	Reduced Mode Downflow Voltage	30
P10	Reduced Mode Inflow Voltage	30
P13	Blower Performance Factor Offset	3/4' = 12 / 5/6' = 10
P14	Blower Performance Factor Window	3/4' = 10 / 5/6' = 5
P15	Silence Alarm Repeat Time (0-60 Minutes)	0
P16	Hold Time for Buttons in Seconds	5
P17	Duration in minutes to exit the service menu	15
P19	Downflow Velocity Increase Display	0
P21	Audible Signal Delay	300
P22	Reset Total Operating Hours	N/A
P23	Audible Alarm Delay for Timer Start	25
P24	Downflow Blower Limit Value Monitoring	50
P25	Current Limit Sash Up	2.5
P26	Current Limit Sash Down	2.5
P27	Auto Alarm Limit	150
P28	Speed Pressure Display	0
P29	Value Window Type 1 Auto Calibration	200
P30	Backup Battery Board	0-No / 1-Yes
P31	Value Window Type 2 Auto Calibration	30
P32	Remote Control/Keyboard Foil Option	0-Rem / 1-Rem &Key
P33	Monitor Alarm	0-Auto / 1-Manual

ERRORS		
E1	Supply Pressure Sensor	Check Sensor Circuit
E2	Exhaust Pressure Sensor	Check Exhaust Sensor
E4	BUS Error	Replace Main Board
E5	RAM Error	
E6	Remote / Memory Error	Check Remote / Replace Board
E8	Memory Error	Replace Main Board
E11	BUS Error	

