

	CO2R_μm l	CO2S_μm l	H2OR_mm l	H2OS_mm l
a	Leaf Fan Speed			15.872
b	current = 2			48.03
	off	1 (Slw)	2 (Fst)	3 4 5
c	0.000257	2.27E-07	-1.32E+03	3.88E-06
	Leaf Fan	* Flow →	Mixer	Temp
	2 Fast	500 μm	OFF	OFF
				Lamp
				OFF