

(B) F l o w C o n t r o l S t a t u s

	F l o w m e t e r	I n p u t P	C h a m b e r P
Type :	s e t p o i n t	s e t p o i n t	s e t p o i n t
S e t P t :	5 0 0 . 0	2 . 0	0 . 0
P c t :	1 0 0 . 0	1 0 0 . 0	1 0 0 . 0
	4 9 9 . 8 u m	2 . 0 1 k P a	0 . 0 2 k P a
P u m p m v :	s t d = 4 5 0 0 ,	m x r = 4 5 0 0 ,	n o w = 4 5 0 0