

Environmental Controls

▼ FLOW: Flow Rate

Control = Flow μm

Target = 500 $\mu\text{mol} / \text{s}$

ΔP to Ref (kPa) = 2

→ ΔP from Chamber (kPa) = 0.02

T) reTarget, exit O) Off, exit



◆ Edit



◆ cancel



Keep