

Environmental Controls

▼ FLOW: Manual

→ { Control = Manual _ %

Target = % F = 100 % I = 100 % C = 100

T) reTarget, exit O) Off, exit

▼ / ► ♦ Edit ♦ ♦ cancel ♦ Keep