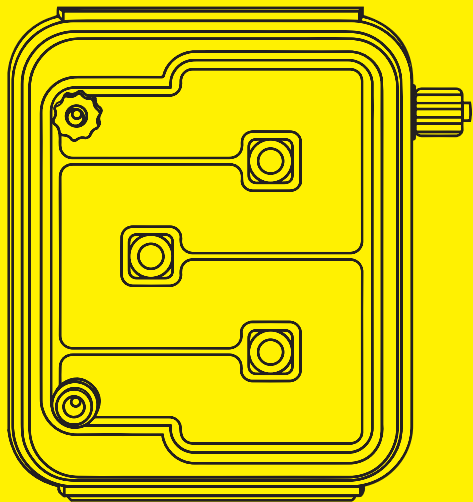


# 4 Ways Vivaflow® SU Will Help You Master TFF Effortlessly



Intuitive setup in **< 5 min**  
ready to process up to 1 L feeds



Runs with a standard pump to **minimize complexity and cost**



Achieves up to 99% recovery **without process optimization**

**Prevents carry-over** while cutting waste by **30.5%**

**SARTORIUS**

Learn More:

