



Clarification Technologies Matrix

WCW

3%

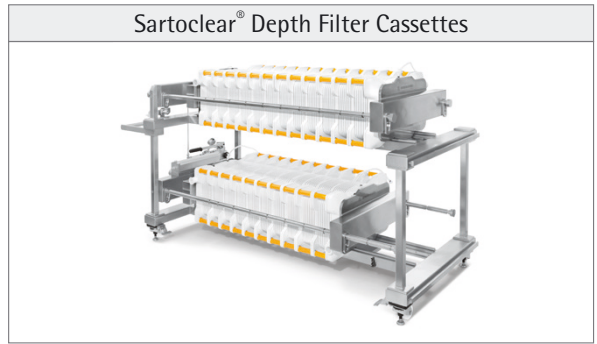
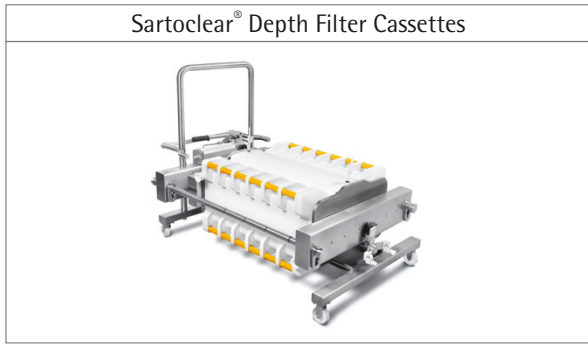
8%

Cell Density ($\times 10^6$ cell/mL)

10

20

Scale
2.000 L
200 L
20 L
2 L



One Step Clarification

Two Step Clarification

Sartoclear® Depth Filters

Sartoclear® Depth Filters

* Volume ranges are dependant on cell & molecule characteristics. Intended as guideline only.



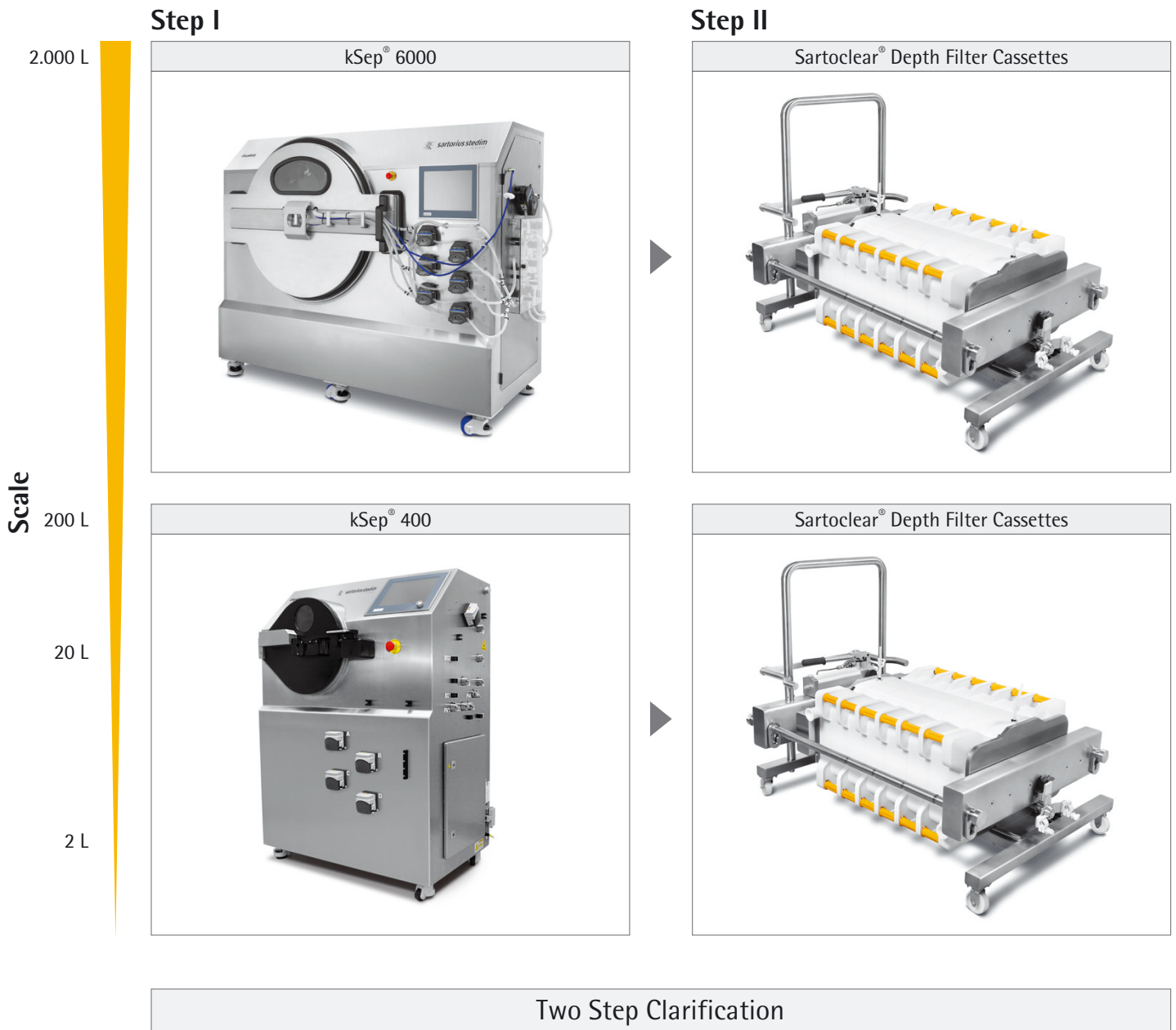
Clarification Technologies Matrix

WCW

> 8%

Cell Density ($\times 10^6$ cell/mL)

> 20



kSep® Centrifugation followed with Sartoclear® Depth Filters

* Volume ranges are dependant on cell & molecule characteristics. Intended as guideline only.