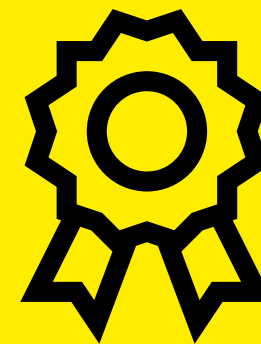
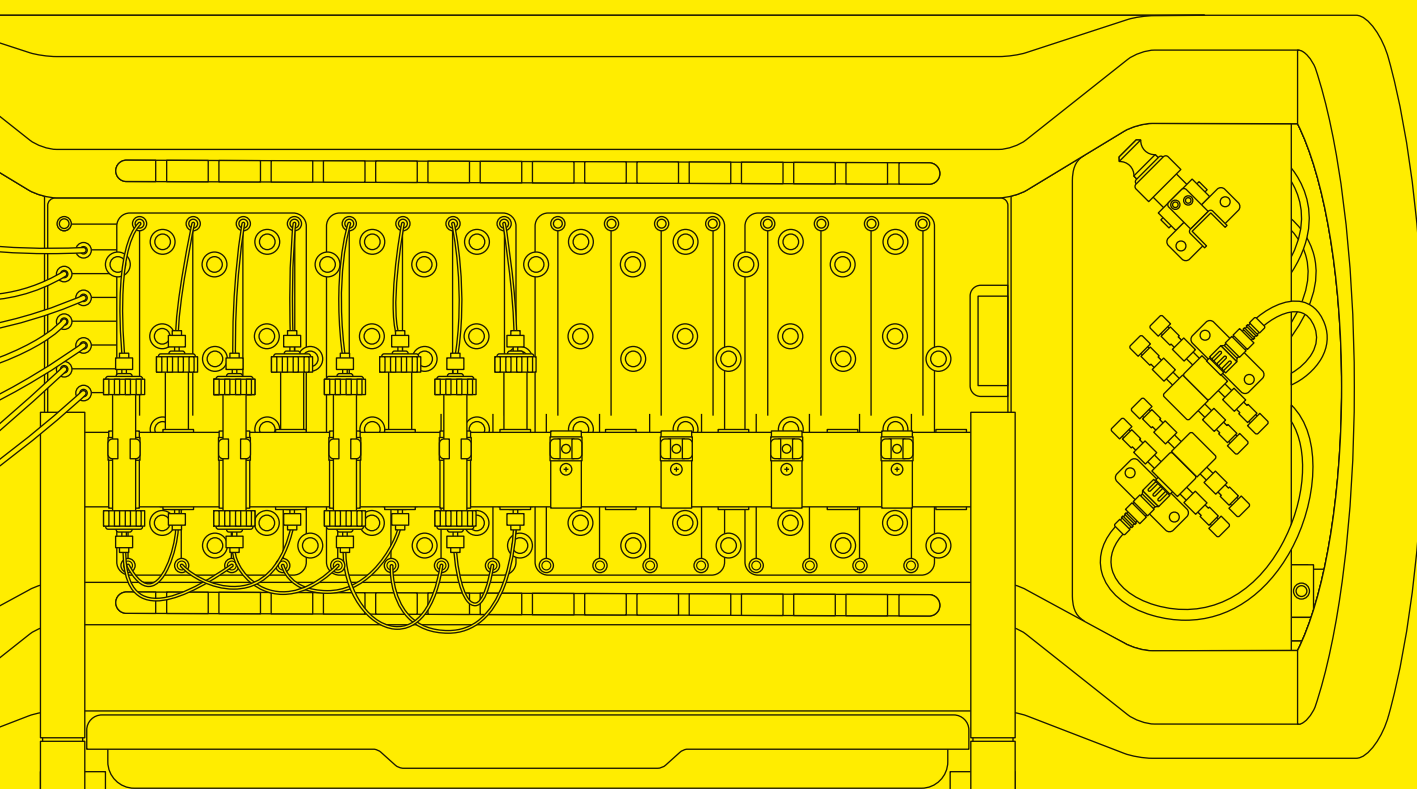


# 5 Reasons to Switch to Multi-Column Chromatography

Up to **80%** savings  
on chromatography resin costs



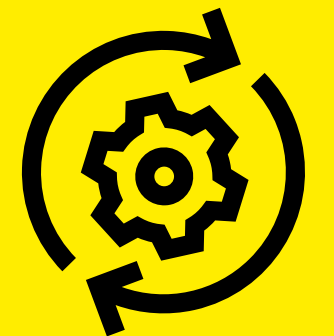
**Critical Quality Attributes  
maintained** when converting  
a batch process



For more information visit:  
[www.sartorius.com/multi-column-chromatography](http://www.sartorius.com/multi-column-chromatography)

Increase productivity  
by **3-5x**

**Easy integration**  
into existing batch  
process



Up to **40%** increase  
in utilization of resin capacity

**SARTORIUS**