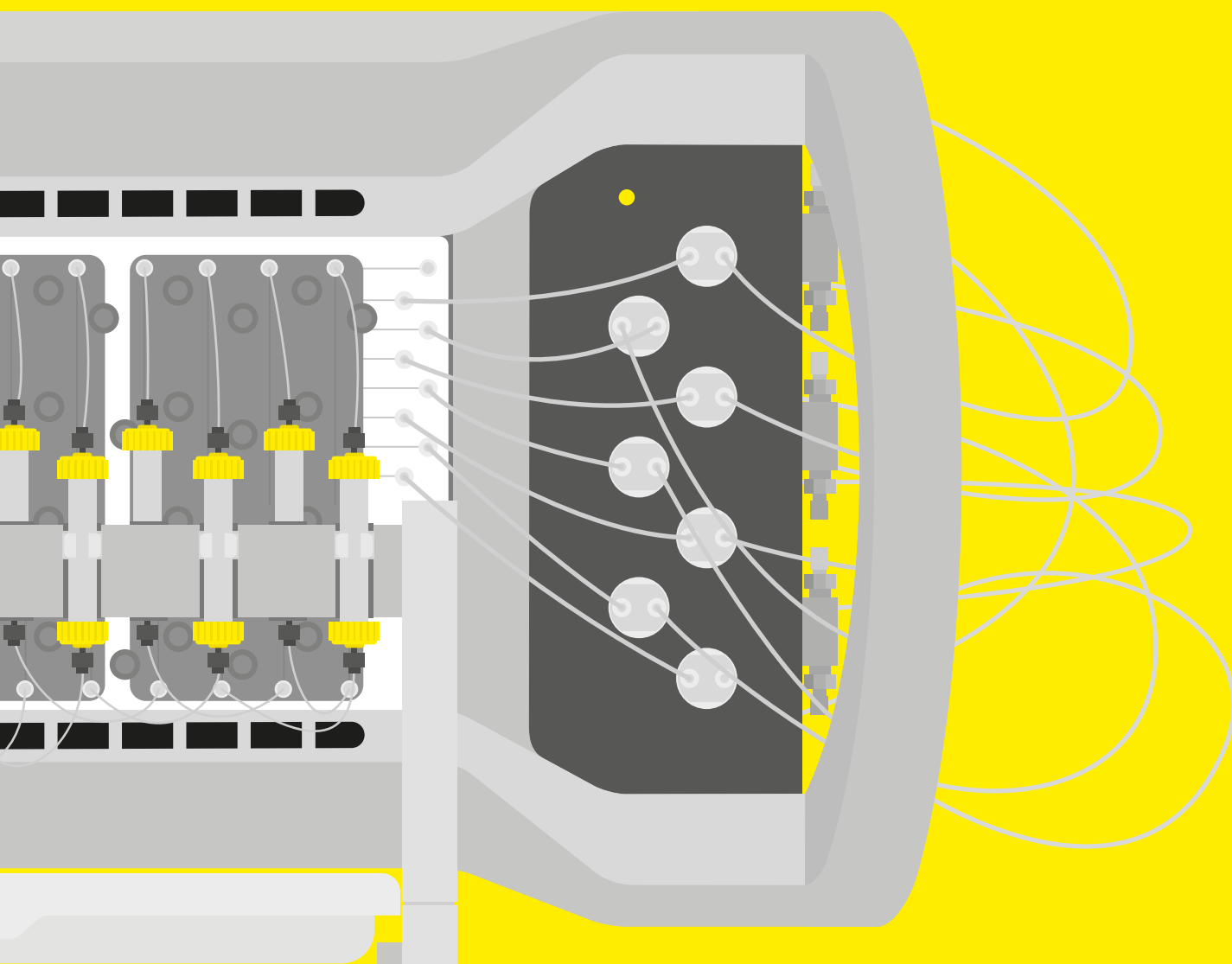
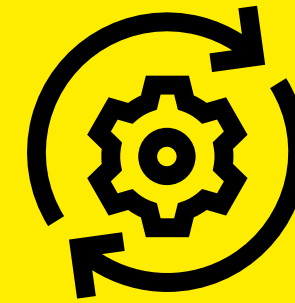


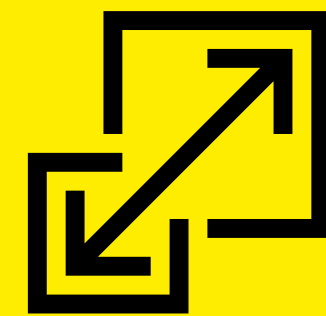
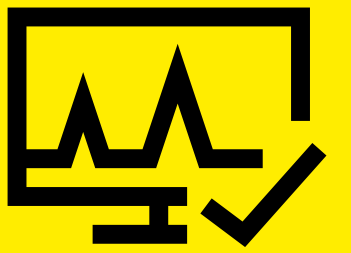
# The Improved Resolute<sup>®</sup> BioSMB PD System



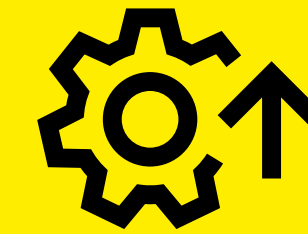
Updated UV-VIS sensor  
technology with  
**maximum robustness**



Same benefits,  
**increased reliability**



**Simplify scalability  
to manufacturing**  
with automation  
platform identical  
to Resolute<sup>®</sup>  
BioSMB 80 | 350

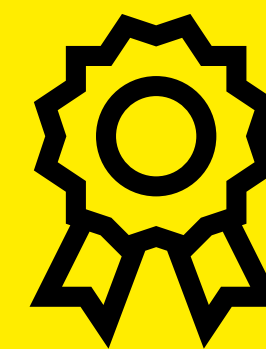


**Increase** productivity  
by 3–5x

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



**Critical quality  
attributes maintained**  
when converting a  
batch process



**Easy integration**  
into existing batch  
process

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

**SARTORIUS**