



## Sartorius PSC Pilot Solution

Data Sheet Collection

**SARTORIUS**