

Success Story: Integrating Vivaspin® as an OEM Component of a POC Test Kit Enables Reliable Diagnostics

Sartorius was recently approached by a third party whose mission is to transform complex patient data into an understandable diagnosis, to inform treatment of metabolic and inflammatory diseases like cancer and obesity. A leader in the field of *in vitro* diagnostics, they make extensive use of Capillary Electrophoresis (CE).

The Challenge

- A new point-of-care (POC) detection kit lacked the sensitivity to detect low levels of the target disease markers
- Disease markers needed to be concentrated without introducing contaminants that could interfere with the diagnostic test
- Reliable supply with global reach was also needed, to support a rapid roll-out of the POC detection kits

The Solution

- Vivaspin® for general laboratory use (GLU)
- High disease marker recovery: PES with suitable MWCO ensures low binding and minimal losses
- Compatible with existing equipment: standard dimensions with conical base works in popular centrifuges
- Easy retentate collection: angular dead stop prevents over-concentration and simplifies sampling for CE

The Value

- Efficient sample concentration increased the sensitivity of CE
- Steady supply secured by regular follow-ups and mid- to long-term forecasts
- Sartorius' filtration expertise allowed the manufacturer to focus on their mission to deliver simple and reliable POC detection kits.



For more information,
visit www.sartorius.com

SARTORIUS