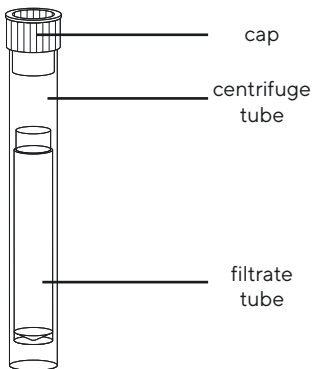


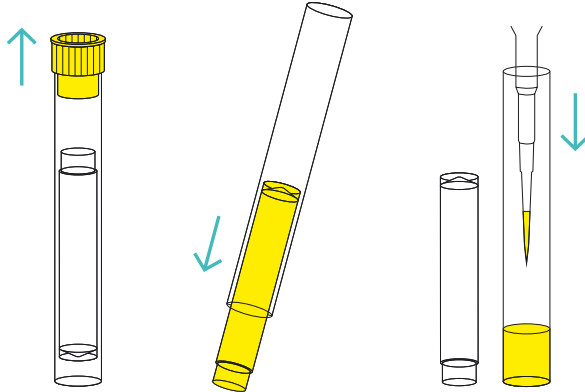
Quick Start Guide Vivaspin® Filtrate

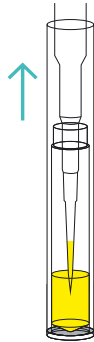
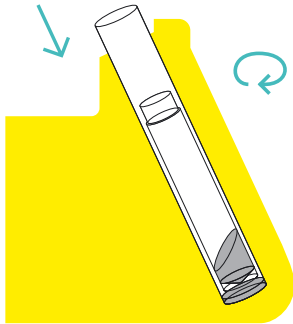
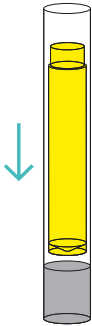


Scan the QR code or search for the order number of your product on www.sartorius.com, to download and read the full Instructions for Use

Vivaspin® Filtrate

- Do **not** use the product after its expiry date.
- Do **not** overfill the filtrate tube.
- Only use the product in suitable centrifuges. Do **not** exceed the RCF limit of the product.







15 mL
17 mm Ø



Centrifuge requirements



Max. filling volume 2.5 mL

Centrifuge with fixed angle rotor $\geq 25^\circ$

RCF limit 2,000 g



Max. filling volume 2.5 mL

Centrifuge with swing bucket rotor

RCF limit 2,500 g
