

Find Your Optimal Centrifugal Concentrator

Just search “Vivaspin”

Centrisart 1 has a new name:
Vivaspin® Filtrate.

By joining the Vivaspin® family, it’s
now easier than ever to find the best
centrifugal concentrator for each of your
ultrafiltration processes.

Now the most comprehensive range
of lab ultrafiltration devices feels more
comprehensive.



Find Your Optimal Centrifugal Concentrator

Vivaspin® Filtrate

Membrane	12 pc
5 kDa CTA	13229-E
10 kDa CTA	13239-E
20 kDa CTA	13249-E
300 kDa PES	13279-E

Find out more:

