

## 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  |
| 100 kDa PES | 13269-GE |
| 300 kDa PES | 13279-E  |



Find out more:

