

Find Your Optimal Centrifugal Concentrator

Just search “Vivaspin”

Vivacell 100 has a new name:
Vivaspin® 100.

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® 100

Membrane	2 pc	10 pc
5 kDa PES	VC1011	VC1012
10 kDa PES	VC1001	VC1002
30 kDa PES	VC1021	VC1022
50 kDa PES	VC1031	VC1032
100 kDa PES	VC1041	VC1042
300 kDa PES	VC1051	VC1052
1,000 kDa PES	VC1061	VC1062
0.2 µm PES	VC1071	VC1072

Find out more:

