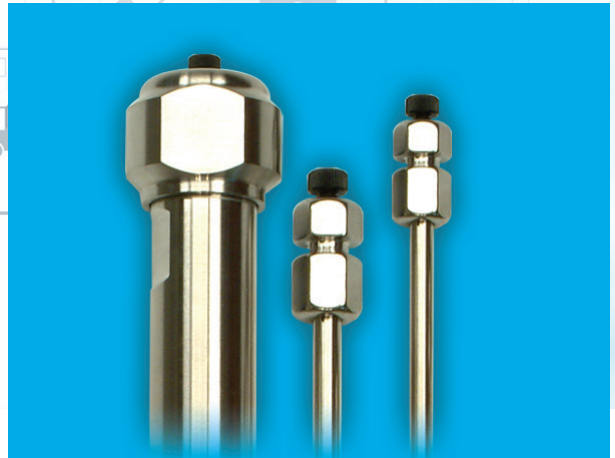


Liquid Chromatography



Spherisep HPLC Columns

Our Spherisep™ line is an equivalent to Waters Spherisorb®. The most popular phase is the ODS2. PerkinElmer offer the Spherisep ODS2 column with equivalent selectivity in 3, 5 or 10 µm particles. Other Spherisep phases are also available including C1, C6, C8, cyano, NH2, ODS1, ODSB, Phenyl, SAX and Silica. Excellent efficiencies, peak shape and resolution are obtained for virtually all Waters Spherisorb HPLC applications from high quality Spherisep HPLC columns. A range of particle sizes offers amazing versatility for capillary and LC/MS to prep and process scale applications.

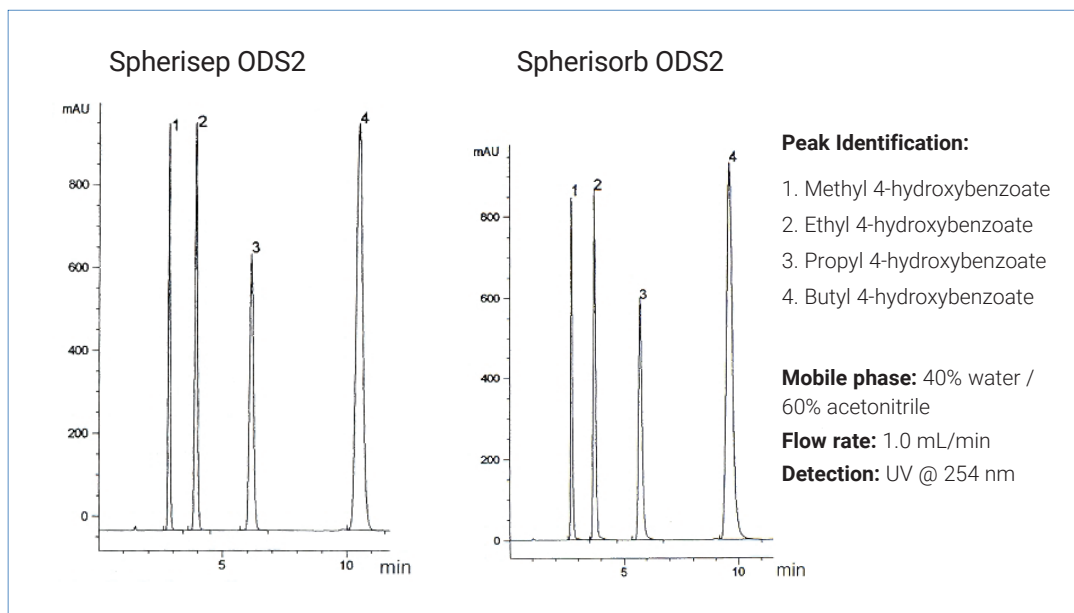


Figure 1: HPLC analysis of 4-hydroxybenzoates using Spherisep ODS2 (left) and Spherisorb ODS2 (right), 150 x 4.6 mm, 5 µm.

Table 1: Spherisep ODS1 HPLC Columns.

Phase	Length (mm)	ID (mm)	Particle Size (µm)	Part No.
Spherisep ODS1	100	4.6	5	125271-SSP-ODS1
Spherisep ODS1	125	4.0	3	104171-SSP-ODS1
Spherisep ODS1	125	4.6	5	105271-SSP-ODS1
Spherisep ODS1	150	4.0	5	134271-SSP-ODS1
Spherisep ODS1	150	4.6	3	135171-SSP-ODS1
Spherisep ODS1	150	4.6	5	135271-SSP-ODS1
Spherisep ODS1	200	4.6	10	145371-SSP-ODS1
Spherisep ODS1	250	3.0	4	153871-SSP-ODS1
Spherisep ODS1	250	4.0	10	154371-SSP-ODS1
Spherisep ODS1	250	4.0	5	154271-SSP-ODS1
Spherisep ODS1	250	4.6	5	155271-SSP-ODS1
Spherisep ODS1	300	4.0	5	164271-SSP-ODS1
Spherisep ODS1 Analytical Guard Cartridges (Pkg. 5)	10	2.0	5	500103-SSP-ODS1
Spherisep ODS1 Analytical Guard Cartridges (Pkg. 5)	10	3.0	5	500101-SSP-ODS1
Analytical Guard Cartridge Holder with integrated coupler	–	–	–	ES500100

Table 2: Spherisep ODS2 HPLC Columns.

Phase	Length (mm)	ID (mm)	Particle Size (µm)	Part No.
Spherisep ODS2	30	4.6	5	185271-SSP-ODS2
Spherisep ODS2	50	4.6	3	115171-SSP-ODS2
Spherisep ODS2	50	4.6	5	115271-SSP-ODS2
Spherisep ODS2	100	2.1	3	122171-SSP-ODS2
Spherisep ODS2	100	4.0	10	124371-SSP-ODS2
Spherisep ODS2	100	4.6	3	125171-SSP-ODS2
Spherisep ODS2	100	4.6	5	125271-SSP-ODS2
Spherisep ODS2	120	4.6	3	1D5171-SSP-ODS2
Spherisep ODS2	125	3.0	3	103171-SSP-ODS2
Spherisep ODS2	250	4.0	5	154271-SSP-ODS1
Spherisep ODS2	250	4.6	5	155271-SSP-ODS1
Spherisep ODS2	125	4.0	5	104271-SSP-ODS2
Spherisep ODS2	125	4.6	3	105171-SSP-ODS2
Spherisep ODS2	150	4.0	3	134171-SSP-ODS2
Spherisep ODS2	150	4.0	5	134271-SSP-ODS2
Spherisep ODS2	150	4.6	3	135171-SSP-ODS2
Spherisep ODS2	150	4.6	5	135271-SSP-ODS2
Spherisep ODS2	250	4.0	5	154271-SSP-ODS2
Spherisep ODS2	250	4.6	10	155371-SSP-ODS2
Spherisep ODS2	250	4.6	3	155171-SSP-ODS2
Spherisep ODS2	250	4.6	5	155271-SSP-ODS2
Spherisep ODS2	300	3.9	5	16e271-SSP-ODS2
Spherisep ODS2 Analytical Guard Cartridges (Pkg. 5)	10	3.0	5	500101-SSP-ODS2
Analytical Guard Cartridge Holder with integrated coupler	–	–	–	ES500100

Other column dimensions, particle sizes, and guard cartridges are available.
Please enquire for more details at LCA.TechSupport@perkinelmer.com