



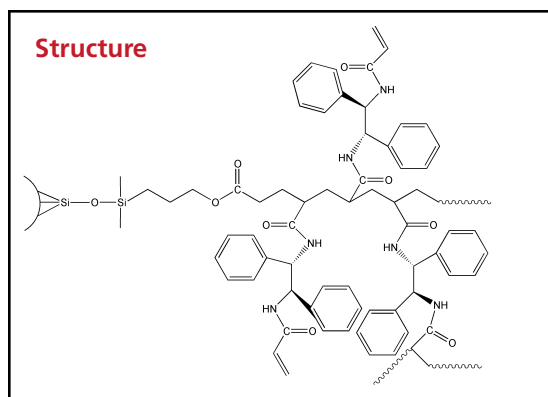
P-CAP-DP™

(R,R) and (S,S)

Poly-CyclicAmine Polymer Chiral Stationary Phase

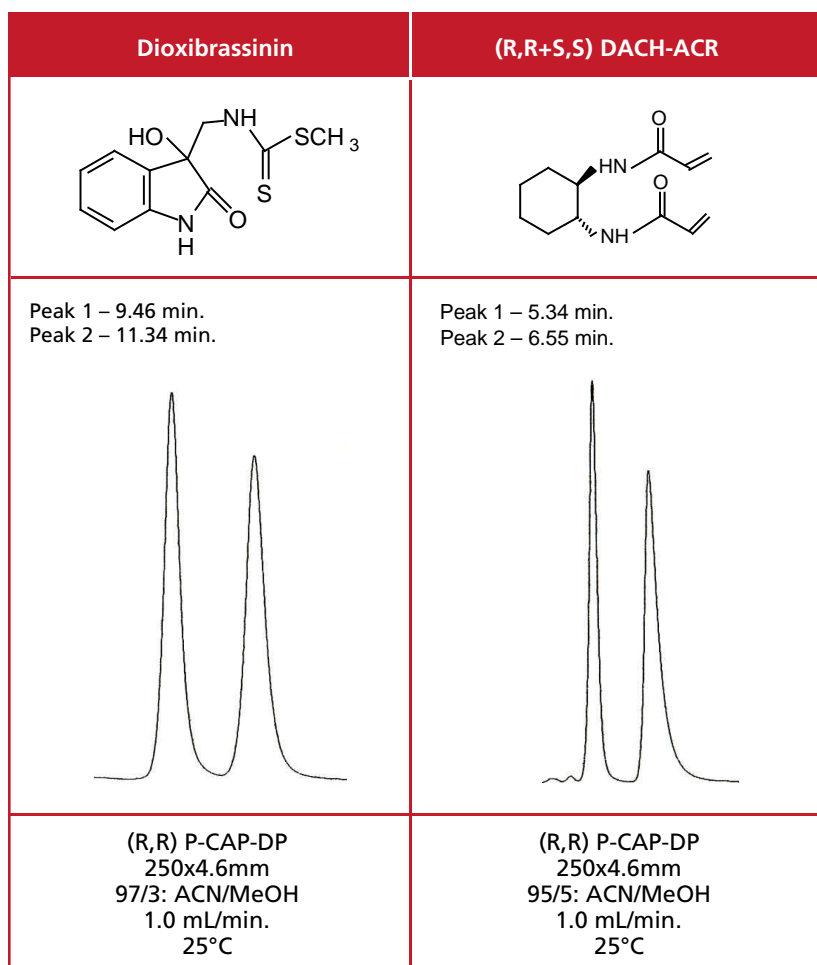
New complementary P-CAP chiral stationary phases with an expanded number of racemates resolved with high selectivity including hydroxycarboxylic acids, alcohols, sulfoxides, esters, amides and lactones and N-blocked amino acids.

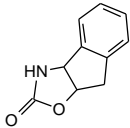
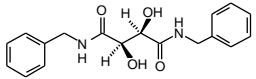
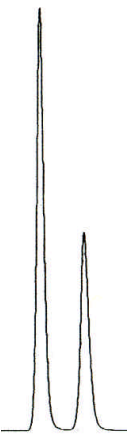
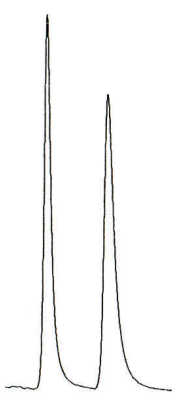
n	Covalently bonded polymer based chiral stationary phase
n	High stability, no memory effect
n	High sample loadability, easy scale-up
n	Similar P-CAP protocols for normal and polar organic phase optimization
n	Normal and polar organic mode chiral stationary phase
n	Reversal of elution order, (R,R) to (S,S) configuration
n	Less polar than P-CAP, ideal for SFC applications



P-CAP-DP expands the capabilities and is complementary to the original P-CAP chiral stationary phase for the normal phase separation of racemic compounds. The introduction of the system, in combination with the chiral hydrogen bonding capabilities offers new opportunities for chiral selectivity. The additional bulk from the aromatic rings further enhances steric effects.

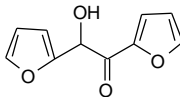
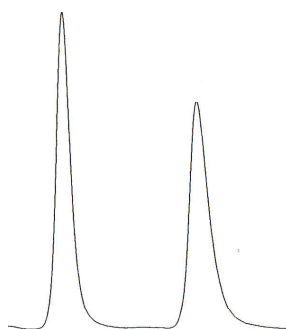
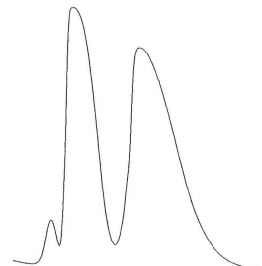
The two basic mobile phase systems employed are the typical heptane/IPA or ethanol and the polar organic mode employing acetonitrile/methanol. Volatile acids and buffers can be used to enhance peak efficiency when needed or to enhance ion detection for MS platforms. There are no known limitations on the kind of solvents that can be used with these phases including halogenated. The well ordered polymer layer is unaffected by solvent changes and allows for excellent efficiency and reproducibility.



<i>cis</i> -(±)-3,3a,8,8a-Tetrahydro-2H-indeno[1,2-d]oxazol-2-one	(±)-N,N'-Dibenzyl-tartaric diamide
	
Peak 1 – 4.87 min. (S) Peak 2 – 5.76 min. (R)	Peak 1 – 5.39 min. (-) Peak 2 – 7.37 min. (+)
	
(R,R) P-CAP-DP 250x4.6mm 98/2: ACN/MeOH 1.0 mL/min., 25°C	(R,R) P-CAP-DP 250x4.6mm 95/5: ACN/MeOH 1.0 mL/min., 25°C

For preparative applications the potential choice of solvents can enhance peak capacity and solubility providing greater sample throughput.

Preparative Application (R,R) P-CAP-DP

Furoin	
	
Load: 8 µg	Load: 1000 µg
Peak 1 – 10.48 min. Peak 2 – 13.20 min.	Peak 1 – 9.45 min. Peak 2 – 11.22 min.
	
(R,R) P-CAP-DP, 250x4.6mm 20/80: EtOH/Hexane 1.0 mL/min. Inj: 0.2 µL (4 µg/µL)	(R,R) P-CAP-DP, 250x4.6mm 20/80: EtOH/Hexane 1.0 mL/min. Inj: 50 µL (20 µg/µL)

AVAILABILITY

Catalog No.	Description	Size
34023	(R,R) P-CAP-DP, 3µm	150x4.6mm
35019	(R,R) P-CAP-DP, 5µm	150x2.1mm
35023	(R,R) P-CAP-DP, 5µm	150x4.6mm
35024	(R,R) P-CAP-DP, 5µm	250x4.6mm
35034	(R,R) P-CAP-DP, 5µm	250x10.0mm
35044	(R,R) P-CAP-DP, 5µm	250x21.2mm
35100	(R,R) P-CAP-DP Guard	2cmx4.0mm
21150	Guard Holder	
36023	(S,S) P-CAP-DP, 3µm	150x4.6mm
37019	(S,S) P-CAP-DP, 5µm	150x2.1mm
37023	(S,S) P-CAP-DP, 5µm	150x4.6mm
37024	(S,S) P-CAP-DP, 5µm	250x4.6mm
37034	(S,S) P-CAP-DP, 5µm	250x10.0mm
37044	(S,S) P-CAP-DP, 5µm	250x21.2mm
37100	(S,S) P-CAP-DP Guard	2cmx4.0mm
21150	Guard Holder	

astec

Advanced Separation Technologies Inc.

World Headquarters: 37 Leslie Court, Post Office Box 297, Whippany, NJ 07981 USA
Tel: 973.428.9080 Fax: 973.428.0152 E-mail: info@astecusa.com www.astecusa.com

UK and Ireland Sales Office: 1 Blake Street, Congleton, Cheshire CW12 4DS UK
Tel: +44 (0) 1260 276276 Fax: +44 (0) 1260 290067 E-mail: info@asteceuro.com