

Table 2. Screening Protocol Leveraging CSPs Selectivity and Mobile Phase Flexibility

Mobile phase system	Mobile phase system success rate	Composition	CSP success rate									Cyclofructans*					
			Polysaccharides			Synthetic			Macrocyclic			Cyclodextrins			Cyclofructans*		
			61%			41%			69%			31%					
			Kromasil CelluCoat	Kromasil AmyCoat	Astec Cellulose DMP	Kromasil TBB	Kromasil DMB	(R,R) P-CAP-DP	Astec CHIROBIOTIC V2	Astec CHIROBIOTIC T	Astec CHIROBIOTIC TAG	Astec CYCLOBOND I 2000	Astec CYCLOBOND I 2000 DMP	Astec CYCLOBOND I 2000 HP-RS	LARIHC CF6-P	LARIHC CF6-RN	LARIHC CF6-M
Normal phase (NP)	44%	NP1	80:20, heptane: isopropanol (with 0.1% TEA and 0.1% TFA)														
		NP2	50:25:25, heptane: isopropanol (with 0.1% TEA and 0.1% TFA):MTBE														
		NP3	80:20:0.1, heptane:ethanol:trifluoroacetic acid														
Polar Organic Mode (POM)	7%	POM1	95:5:0.3:0.2, acetonitrile:methanol: acetic acid:triethylamine														
		POM2	95:5, acetonitrile:isopropanol (with 0.1% TEA and 0.1% TFA)														
		POM3	60:40:0.3:0.2, acetonitrile:methanol:acetic acid:triethylamine														
Polar Ionic Mode (PIM)	17%	PIM	100:0.1:0.1, methanol:acetic acid:triethylamine														
Reversed-Phase (RP)	28%	RP	50:50, 20 mM ammonium acetate (pH 4.0):methanol														

* The cyclofructan phases have found most success with chiral primary amines and are therefore screened only on those compounds.