Polymer & Solvent Typical samples Solvent Analysis temp Column(s) Standards Bottles, bags, pipes, car fenders, Polyolefins, polyethylene, Trichlorobenzene 140 to 180 °C PLgel Olexis Polystyrene PLgel MIXED-A polypropylene PLgel MIXED-B Water (pH adjusted), buffers Ambient to 40 °C PL aquagel-OH lonic polymers, chitosan, Polyethylene glycol/oxide insecticides, slow release agents, polyethylene glycol, polyacrylic acid, polyvinyl pyrrolidone DMS0 + 0.1% LiBr 50 to 80 °C PLgel MIXED-A Vater treatment, agrochemical Lignin Polyethylene glycol/oxide PLgel MIXED-B Thickeners and stabilizers DMS0 + 0.1% LiBr 50 to 80 °C PLgel MIXED-A Polyethylene glycol/oxide PLgel MIXED-B hickeners, slow release energy Low MW cellulose, Water (pH adjusted), buffers Ambient to 40 °C PL aquagel-OH Polyethylene glycol/oxide agents, gelling agents, baking aids polysaccharides, dextran, pectin, ackaging, plastic wrap, plastic Vinylic polymers and copolymers, Tetrahydrofuran Ambient to 40 °C PLgel MIXED-C Polystyrene, PLgel MIXED-D polyethylene terephthalate*, polymethylmethacrylate oottles and caps polyvinylidene chloride PLgel MIXED-E Paints, adhesives, binders, art Ambient to 40 °C PLgel MIXED-C Polystyrene, Alkyd, epoxy, acrylic polymers, PLgel MIXED-D polymethylmethacrylate PLgel MIXED-E Phenol and urea formaldehyde, DMSO + 0.1% LiBr, DMF or 50 to 80 °C PLgel MIXED-B Polymethylmethacrylate or Adhesives, molded products, PLgel MIXED-C electrical and heat insulators melamine, cellulose derivatives polyethylene glycol/oxide Fibers, toothbrush bristles, tubing, Polyesters, polyamides, nylons Hexafluoroisopropanol PL HFIPgel Polymethylmethacrylate fishing line, low strength machine aceutical (drug delivery) Hydrophilic controlled, targetted, Polyvinyl pyrrolidone, polylactic Water (pH adjusted), buffers Ambient to 40 °C PL aquagel-OH Polyethylene glycol/oxide acid, polyglycolic acid, polystyrene localized drug release Hydrophobic controlled, targeted, Polydimethyl siloxane, PLgel MIXED-C Polystyrene, localized drug release polyurethane, ethylene vinyl PLgel MIXED-D polymethylmethacrylate acetate PLgel MIXED-E Polyethylene**, polyurethane, Tetrahydrofuran Ambient to 40 °C PLgel MIXED-C Artificial joints, contact lenses, Polystyrene, PLgel MIXED-D polymethylmethacrylate silicones, polylactide/glycolide and copolymers Plumbing pipes, guttering, plastic Polyvinyl chloride, polycarbonate, Tetrahydrofuran Ambient to 40 °C PLgel MIXED-B Polystyrene, PLgel MIXED-C tableware, compact discs, glazing, ABS, polymethylmethacrylate, polymethylmethacrylate PLgel MIXED-D automotive applications Tires, fuel hoses, insulating PLgel MIXED-B Natural and synthetic rubber, Ambient to 80 °C Polystyrene materials, lubricating agents, heat polydimethyl siloxane, PLgel MIXED-C resistant tiles polyethylene wax Special conditions required for each - contact Agilent technical support Polyphenylene sulfide, PTFE, Non-stick coatings, engineering polymers, detergents, PEEK, liquid crystal polymers, www.agilent.com/chem/techsupport high strength and chemical esistant applications

* Refer to "Synthetic fibers" for analysis, ** Refer to "Petrochemical" for analysis

Troubleshooting Solution Minimize tubing length Tighten injection seal Check connector fitting Replace or repair column Column degradation Use mobile phase additives Shear degradation Check age of standards Modify dissolution process (no excessive shaking) Peak broadening Minimize tubing/check fittings Eluent too viscous Heat column oven If possible use smaller cell volume Column failure Replace or repair column Add higher pore size column Baseline drift/noise mn/detector contamination Flush column/detector to waste Clean eluent Use better quality solvents **Bubbles** in detector De-gas solvent Use column heater/insulate tubing Wait for previous run to finish Ensure injection valve is flushed clean Reverse signal polarity of RI dn/dc of polymer less than solvent Change eluent to get true peak Cannot be avoided, but reduced by preparing sample in mobile phase Split peaks Reduce loading/loop size ple loading too large Blocked/partially blocked frit Replace frit - use 2 µm inline filter to stop clogging Replace column Partially blocked injection valve Replace rotor seal Check for bubbles in pump head De-gas solvent Use modifiers/additives Interaction with packing Change eluent polarity High pressure umn/frit blockage Replace frit - use 2 µm inline filter to stop clogging Filter samples Use guard column Reverse flow on column to clear blockage Detector blockage Flush with solvent Heat columns, ensure solvent bottles are not cold Filter samples Pressure fluctuation ild-up of particulates Check valve is dirty Replace/clean check valve Purge pump/de-gas solvent

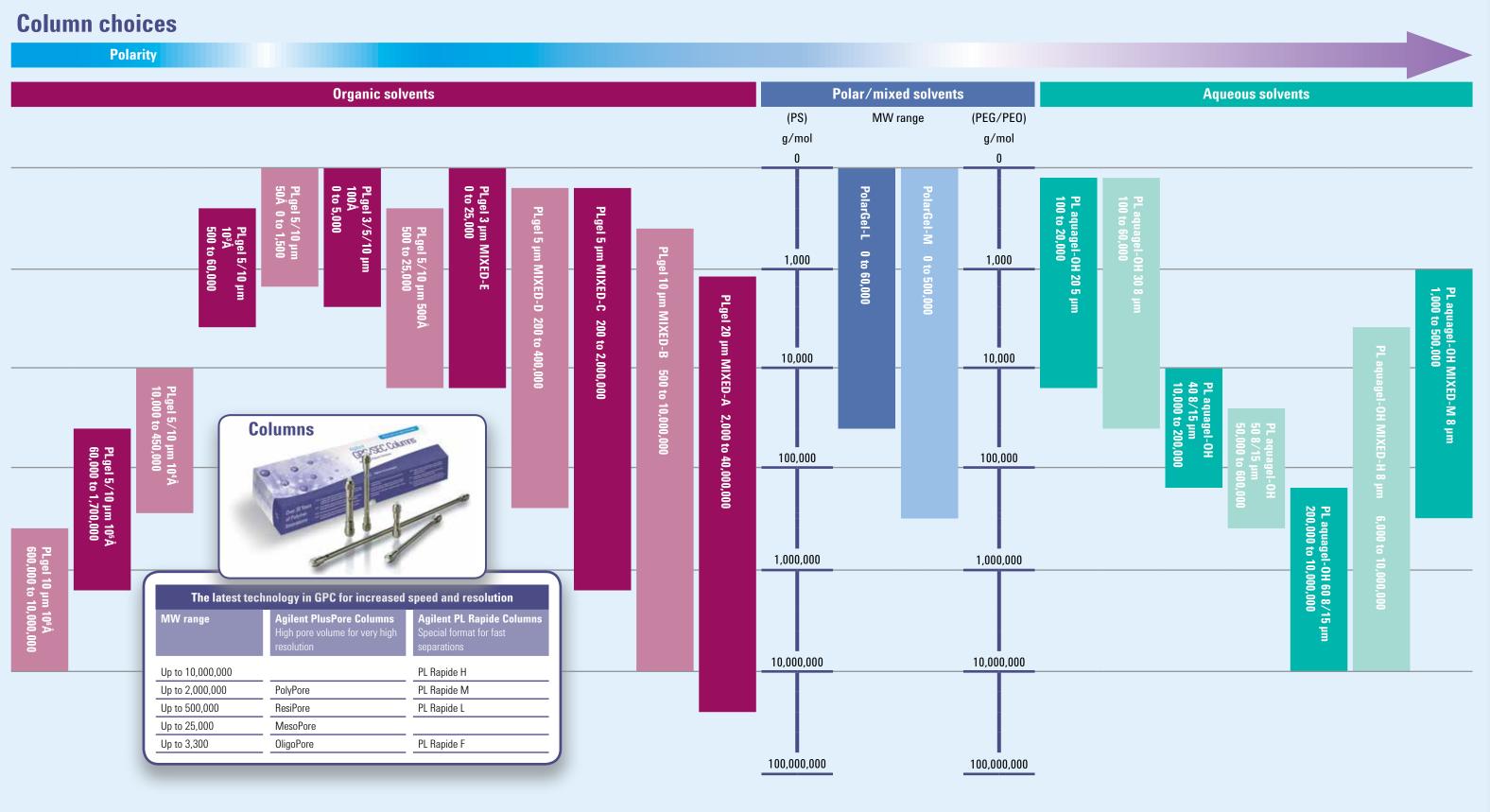
Insufficient flow to pump

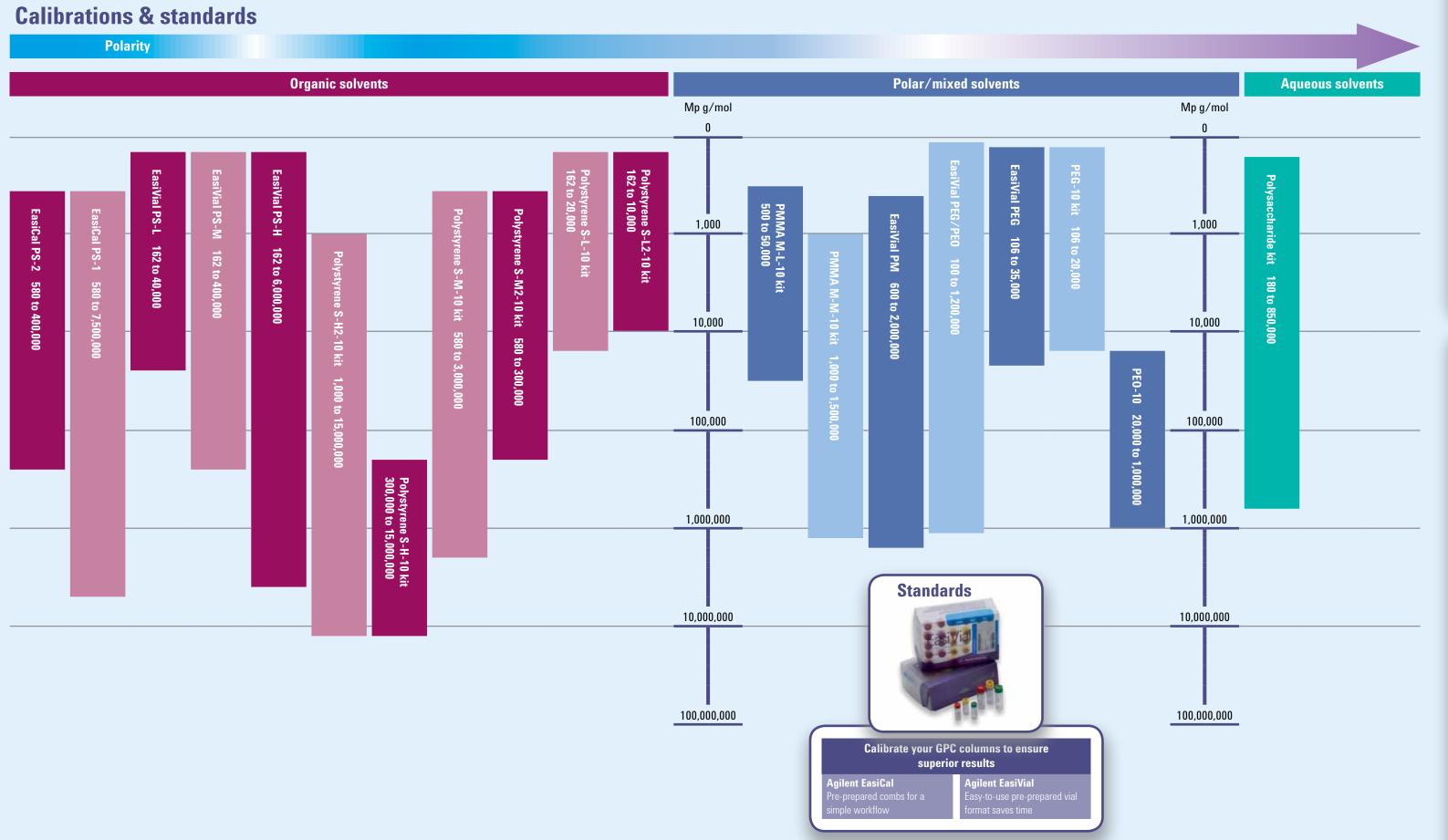
Clean mobile phase inlet

Elevate reservoir above pump head

ACHIEVE MORE WITH THE POLYMER ANALYSIS PEOPLE

The Measure of Confidence





Transfer guide

Organic columns

Flush with new eluent at 0.5 mL/min

30 to 40 °C* as required for analysis

for two column volumes

at 1 °C/min

••••

Column supplied in ethylbenzene

Flush column direct with new eluent**

100 to 220 °C* as required for

Flush with new eluent at 0.2 mL/min

Increase column temperature to

50 to 80 °C* as required for analysis

for two column volumes

required flow rate

Flush column with acetone at Flush column with acetone at Set column oven to 50 °C, flow at 0.5 mL/min for two column volumes 0.5 mL/min for two column volumes

Flush with new eluent at 0.2 mL/min

50 to 80 °C* as required for analysis

at 50 °C at 0.1 mL/min for two for two column volumes column volumes Increase column temperature to Increase column temperature to Increase column temperature to

at 1 °C/min analysis at 1 °C/min Operate column in new eluent at Operate column in new eluent at Operate column in new eluent at

required flow rate required flow rate *Always ensure operating temperature is at least 10 °C below boiling point of solvent.

**Always ensure miscibility. If unsure, use acetone at room temperature.

Polar columns			
Column supplied in water containing 0.02% NaN ₃			
Transfer to aqueous e.g. Water Buffer	Transfer to mixed solvent systems e.g. Water/THF* Water/methanol*	Transfer to polar organic e.g. DMF DMS0	
↓	\	_	
Flush column with pure water at 1.0 mL/min for two column volumes	Flush column with pure water at 1.0 mL/min for two column volumes	Flush column with pure water at 1.0 mL/min for two column volumes	
↓	\downarrow	↓	
Flush with new buffer at 1.0 mL/min for two column volumes	Flush with new, pre-mixed eluent at 0.2 mL/min for two column volumes	Flush with acetone at 0.5 mL/min fo two column volumes	

Increase column temperature* as

Operate column in new eluent at Operate column in new eluent at required flow rate required flow rate

50 °C maximum, using required flow required for analysis at 1 °C/min

at 1 °C/min *Always ensure operating temperature is at least 10 °C below boiling point of solvent.

> Operate column in new eluent at required flow rate

Operate column in new eluent at

Aqueous columns			
Column supplied in water containing 0.02% NaN ₃			
Transfer to aqueous e.g. Water Buffer	Transfer to mixed solvent systems e.g. Water/methanol*		
↓ Flush column with pure water at 1.0 mL/min for two column volumes	↓ Flush column with pure water at 1.0 mL/min for two column volumes		
\downarrow	\downarrow		

Flush with new buffer at 1.0 mL/min Flush with new, pre-mixed eluent at for two column volumes 0.2 mL/min for two column volumes Operate column in new eluent at

Increase column temperature** as 50 °C maximum, using required flow required for analysis at 1 °C/min

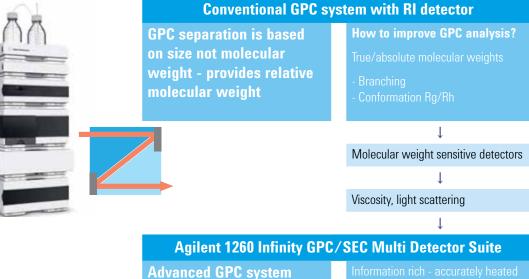
Operate column in new eluent at

Operate column in new eluent at required flow rate

required flow rate *Maximum of 50% organic solvent.

**Always ensure operating temperature is at least 10 °C below boiling point of solvent.

Advanced GPC









Agilent Technologies Published in the UK Feb 27, 2014 www.agilent.com/chem/GPCresources Publication Number: 5991-3802EN