

SELECTING YOUR HPLC COLUMN

The Right Column Makes All the Difference



Low MW
(>5000 g/mol)



Water Soluble

- Ion Exchange
- Reversed Phase
- HILIC



Organic Soluble

- Normal Phase
- Reversed Phase
- Non-Aqueous Size Exclusion



High MW
(<5000 g/mol)



Water Soluble

- Aqueous Size Exclusion
- Ion Exchange
- Hydrophobic Interaction



Organic Soluble

- Non-Aqueous Size Exclusion