







Gas Chromatography Inlets

	Inlet	Column	Mode	Sample Concentration	Sample to Column	GC Compatibility	Comments
 S/SL	Split/Splitless	Capillary	Split Pulsed split Splitless Pulsed splitless	High High Low Low	Very little Very little All All	Intuvo 9000 8890 8860	<ul style="list-style-type: none"> • Most commonly used inlet • Very flexible
 MMI	Multimode Inlet	Capillary	Split Pulsed split Splitless Pulsed splitless Solvent vent Direct	High High Low Low Low	Very little Very little All All Most All	Intuvo 9000 8890	<ul style="list-style-type: none"> • Flexibility of standard S/SL and PTV inlets • Large-volume injections • Cool injections • Controlled speed injections, concentrate analytes, and vent solvent
 COC	Cool On-Column	Capillary	N/A	Low or labile	All	8890 8860	<ul style="list-style-type: none"> • Minimal decomposition and discrimination
 PP	Purged Packed	Packed Megabore capillary	N/A	Any	All	8890 8860	<ul style="list-style-type: none"> • Typically used when high-efficiency separations are not required
 PTV	Programmable Temperature Vaporizing	Capillary	Split Pulsed split Splitless Pulsed splitless Solvent vent	High High Low Low Low	Very little Very little All All Most	8890	<ul style="list-style-type: none"> • Controlled speed injections, concentrate analytes, and vent solvent
 VI	Volatiles Interface	Capillary	Direct Split Splitless	Low High Low	All Very little All	8890	<ul style="list-style-type: none"> • Lowest dead volume • Maximum flow of 100 mL/min • Cannot perform manual injections • Often used for headspace, purge and trap (P&T), and thermal desorption