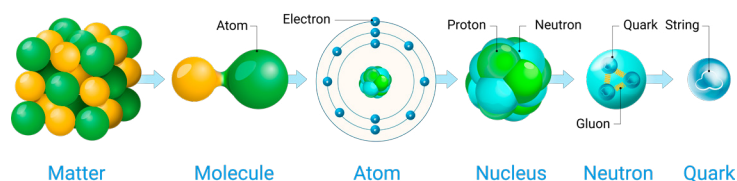


Vacuum for Particle and Plasma Physics



Why you need vacuum	To avoid plasma and particle beam being disrupted by interaction with ambient gases
Typical vacuum level	From 10^{-8} and 10^{-11} mbar
Typical gases	Light gases as He and H ₂
Suitable vacuum technologies	Ion pumps, titanium sublimation combination pumps, turbomolecular pumps, scroll pump, leak detectors
Products	<ul style="list-style-type: none"> – Turbomolecular pumps and turbo pumping systems: TwisTorr 84FS, TwisTorr 305FS, TwisTorr 704FS, Turbo-V 2300, TPS-mobile, TPS-flexy. – Ion and getter pumps: Vaclon Plus Pumps from 150 to 1000 L/s, Titanium Sublimation Combination Ion Pumps (TSP), Non-Evaporable Getter (NEG) Combination Ion Pumps. – Leak detection: Helium Mass Spectrometers HLD. – Vacuum measurement: Inverted Magnetron & Bayard Alpert Ion Gauge tubes.
Vacuum solutions	Multiple pumping stations along the vacuum chamber.
Typical requirements	<ul style="list-style-type: none"> – High Pumping speed for light gases – Pump remote operation (up to 100 m) – Long lifetime – Resistance to mag-field up to 100 Gauss (10 mT) for turbomolecular pumps – Resistance to radiation up to 5×10^5 Gray for turbomolecular pumps