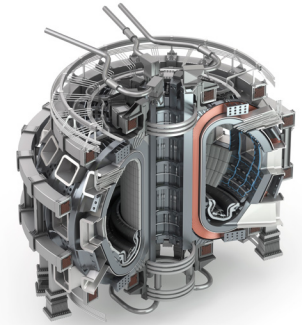


Vacuum for Nuclear Fusion Technology



Why you need vacuum	To create the right environment for atoms fusion, requiring a confined overheated and ultra-clean plasma.
Typical vacuum level	From 10^{-8} and 10^{-11} mbar
Typical gases	Hydrogen, tritium and deuterium
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 100m) - 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