

AGILENT TWISTORR 304 FS

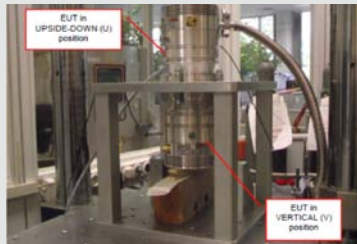
HIGH PERFORMANCE, INNOVATION, RELIABILITY.

Agilent commitment to technology leadership translates into quality, reliability and performance. Our products need to comply with strict parameters, to make sure we can meet and exceed customer expectations. The TwisTorr 304 FS represents a breakthrough in turbomolecular pump technology, based on two patented innovations, TwisTorr stages and Floating Suspension, which strongly impact performance and reliability. Development of the 304 FS included a comprehensive series of tests and analysis, performed to ensure the highest levels of pump performance and reliability. Through those tests VPD verified, step by step, that pump performance and robustness is ready to meet our customers' demanding specifications.

This rigorous test regime of product performance, reliability and durability according to Agilent standards ensures the release of a robust, high performance, system compatible product.



TwisTorr 304 FS Test Elements



Shock Test



Vibration Test



Temperature Test



Noise Test



Reliability assessment on Agilent system



Life Tests



Vacuum Test



Humidity Test



Agilent Technologies

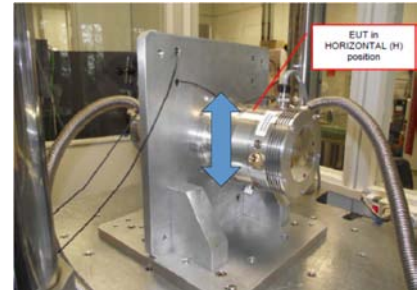
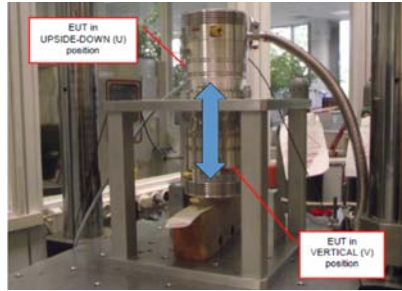
Shock Test

4 Pumps tested in each position.

Operative: Pump tested under nominal working conditions (pump ON). After the test the pump must operate as new.

Non Operative: Pump not running (pump OFF) to simulate transportation and storage. The pump after the test must operate as new.

Pump position	Sinusoidal (Operative)	Trapezoidal (Non-Operative)
Horizontal	> 110g	> 40g
Vertical	> 150g	> 60g
Upside-down	> 150g	> 60g



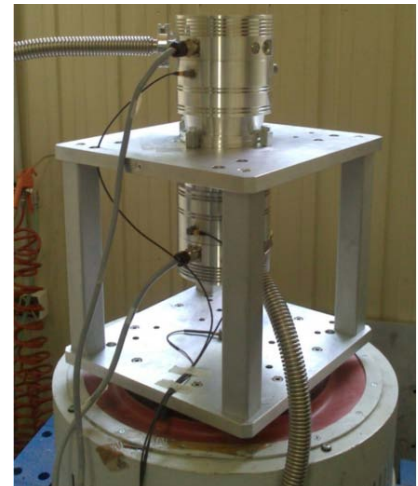
Vibration Test

4 Pumps tested in horizontal, vertical and upside position.

Operative: Pump tested under nominal working conditions (pump ON). After the test the pump must operate as new.

Non Operative: Pump not running (pump OFF) to simulate transportation and storage. The pump after the test must operate as new.

Pump position	Operative (5-500Hz)	Non-Operative (5-500Hz)
Horizontal	0.2grms	2grms
Vertical	0.2grms	2grms
Upside-down	0.2grms	2grms



Temperature Test

4 Pumps tested in horizontal, vertical and upside position.

Operative: Pump tested under nominal working conditions (pump ON). After the test the pump must operate as new.

Non Operative: Pump not running (pump OFF) to simulate transportation and storage. The pump after the test must operate as new.

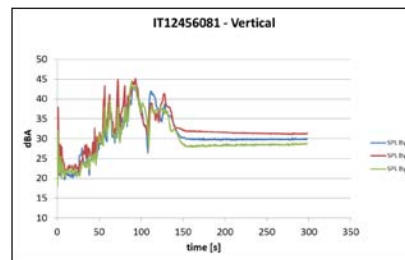
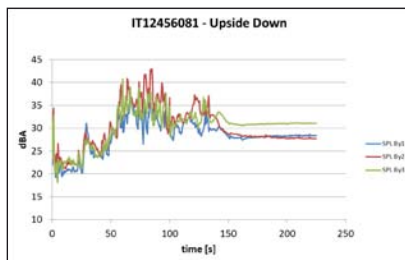
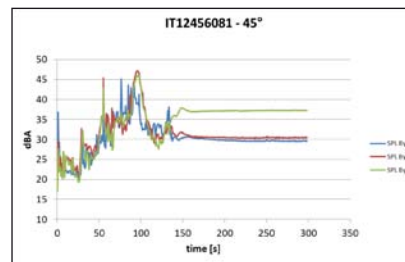
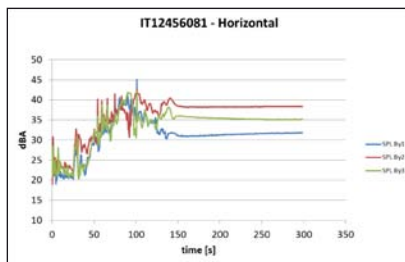
Pump position	Operative	Non-Operative
Horizontal	0 °C to 40 °C	-40 °C to +70 °C
Vertical	0 °C to 40 °C	-40 °C to +70 °C

Entire thermal cycle = 80 hours



Noise Test

Noise specs are measured at full speed at 1 meter from the pump (noise specs < 50dB (A)).

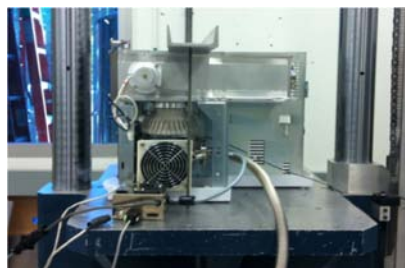


Reliability assessment on Agilent GCMS

Non-operational shock test: system drop with pump not running, after each drop pump is switched on and

after stabilization unbalance, Fourier spectra, power adsorption and bearing temperature are acquired.

# shock	1	2	3	4
TV301 IT12186116	pass	pass	pass	pass
TV301 IT12196067	pass	pass	pass	pass
TV301 IT12166229	pass	pass	pass	fail
304 - LP 13	pass	pass	pass	pass
304 - LP 15	pass	pass	pass	pass
304 - LP 16	pass	pass	pass	pass
304 - LP 14	pass	pass	pass	pass



Operational shock test: system drop with pump running, after each drop unbalance, Fourier spectra, power adsorption and bearing temperature are acquired.



# shock	Increasing drop height																				
	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	...	25	
Height (inches)	3	3	3	3	3	3	3	3	3	3	3.1	3.2	3.3	3.4	3.6	3.8	4	4.3	...	10	
TV301 IT12186116	crash																				
TV301 IT12196067	crash																				
TV301 IT12166229	crash																				
304 - LP 13	pass	pass	pass	pass	crash																
304 - LP 15	pass	crash																			
304 - LP 16	pass	Pump perfectly working but unbalance above baseline																			
304 - LP 14	pass	Pump perfectly working but unbalance above baseline																crash			

Pass = Unbalance within baseline value

