

Oil and gas applications – refinery plant

Vibration sensors improve the reliability and efficiency of oil and gas machines, so you can minimize unplanned downtime, reduce operating costs and protect critical machinery from damage or failure.



Vibration monitoring for oil and gas machinery and processing equipment

Oil and gas operations top the list of extreme operating environments. Detecting and preventing unexpected machinery and component problems is a priority. Vibration monitoring detects baseline and trend machinery health, provides a cost-efficient, reliable choice for maintaining machine health, extending equipment life and maximizing operational efficiency.

Vibration monitoring benefits

- Acceleration and velocity transmitters provide true peak detection
- Low cost implementation
- Increased safety and asset management
- Better alarm handling and reporting
- Programmable alarm functions for automatic work order generation
- Optimized maintenance scheduling
- Hazardous area certified models

Quickly identify machinery faults

- Bearing damage*
 - Lubrication issues*
 - Imbalance*
 - Misalignment
 - Bent shaft
- * primary causes of equipment failure



Typical oil and gas machinery using vibration monitoring solutions

- Blowers
- Centrifuges
- Chillers
- Compressors
- Cooling towers
- Fans
- Gearboxes
- Mixers
- Motors
- Pumps
- Reciprocating compressors
- Turbines
- Vibrating conveyors



Authorised distributor

In Australia:

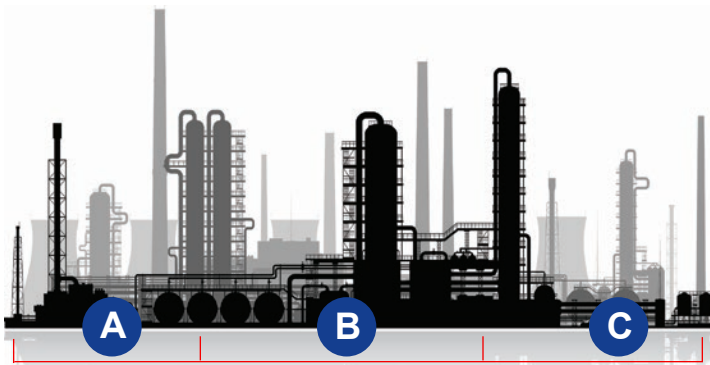
For customer service, call 1300-735-292
To email an order, ordersau@thermofisher.com
To order online: thermofisher.com

In New Zealand:

For customer service, call 0800-933-966
To email an order, ordersnz@thermofisher.com
To order online: thermofisher.com

Reliable vibration monitoring helps you understand asset condition

Sensors for data driven decision-making | Critical vibration measurement points | 1000's of sensors per plant



A Hazardous area
Class I, Div 1
Class I, Zone 0/1

B Explosion-proof

C Class I, Div 2
Class I, Zone 2

Certified hazardous approvals



A = 786A-IS, 787A-IS



B = PC420VP-10-EX
PC420VP-20-EX



C = 786-500-D2, 786A-D2



Key considerations, recommended solutions and products for oil and gas applications

Consideration	Solution	Product
A - Hazardous environment, intrinsically safe	Class 1 Division 1 (Zone 0/1) certified sensors with safety barriers	786A-IS 787A-IS
B - Hazardous environment, explosion-proof	Explosion-proof sensors, fully self-contained	PC420VP-10-EX PC420VP-20-EX
C - Hazardous environment, Class 1, Div 2, Zone 2 Trendable sensors Standard sensors	4-20 mA output sensors Class 1 Division 2, Zone 2 sensors and cables	786-500-D2 786A-D2 786T-D2 R6D2-0-J9T2A/J9T3A

Why choose Wilcoxon?

- Customer support
- Reliability and quality
- Consistent, long-term product performance
- One-stop for everything you need

Complete solutions to meet your vibration monitoring needs!

Specialty and general purpose sensors, cables and connectors, mounting accessories, enclosures, portable meters, intelligent transmitters

Authorised distributor

In Australia:

For customer service, call 1300-735-292
To email an order, ordersau@thermofisher.com
To order online: thermofisher.com

In New Zealand:

For customer service, call 0800-933-966
To email an order, ordersnz@thermofisher.com
To order online: thermofisher.com