

Pump monitoring guide



Key indicators of pump health

- Balance
- Alignment
- Cavitation
- Vane pass frequency
- Bearing health



4 common reasons pumps fail

- Impeller wear
- Insufficient lubrication
- Improper sizing
- Seal wear



Protect your pumps

- Identify faults early
- Pump cavitation - detect and correct process
- Extend equipment life
- Cost-effective predictive maintenance
- Minimize lost production time
- Ensure maximum efficiency of plant operations



Pumps are crucial components of many processes across most industrial applications. Vibration monitoring is a cost-effective way to maintain pump and overall machine health, minimize downtime and reduce operating costs.

Recommended monitoring solutions

Consideration	Solution	Product
Pump cavitation	High frequency sensor	786A PC420-ATP iT300
Underwater pumps	Watertight sensor	786F 786/R6Q connector
Contamination, misalignment, bearing failure	Dynamic vibration monitoring	786A 786A and iT300

Ideal markets and applications

- Oil and gas
- Power generation
- Pulp and paper mills
- Steel mills
- Food processing
- Plastic manufacturing

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