

Cement manufacturing applications

Vibration sensors help improve reliability and efficiency so you can minimize unplanned downtime, reduce operating costs and protect critical components from damage or failure.



Vibration monitoring for cement manufacturing

Using accelerometers and 4-20 mA sensors for condition monitoring in cement manufacturing plants is proven to maximize operational uptime and minimize maintenance costs. Detecting and preventing unexpected machine failures is a priority. Vibration monitoring helps establish machine baselines, allowing trend histories of key components. This provides a cost-saving, reliable choice for maintaining components, extending equipment life and maximizing operational efficiency.



Vibration monitoring benefits

- Vibration levels can reveal machine's health and offer insight into developing faults
- Low-cost implementation – can be easily incorporated into existing control systems
- Increased safety and asset management
- Better alarm handling and reporting
- Programmable alarm functions for automatic work order generation
- Optimized maintenance scheduling

Monitor vibration levels on critical pieces of equipment

- Motor bearings
- Air compressors
- Belt drives
- Kiln rollers
- Fans and blowers
- Hammer mills



Quickly identify machinery faults

- Bearing damage*
- Misalignment*
- Imbalance*
- Contamination
- Bent shaft

* primary causes of equipment failure

Authorised distributor

In Australia:

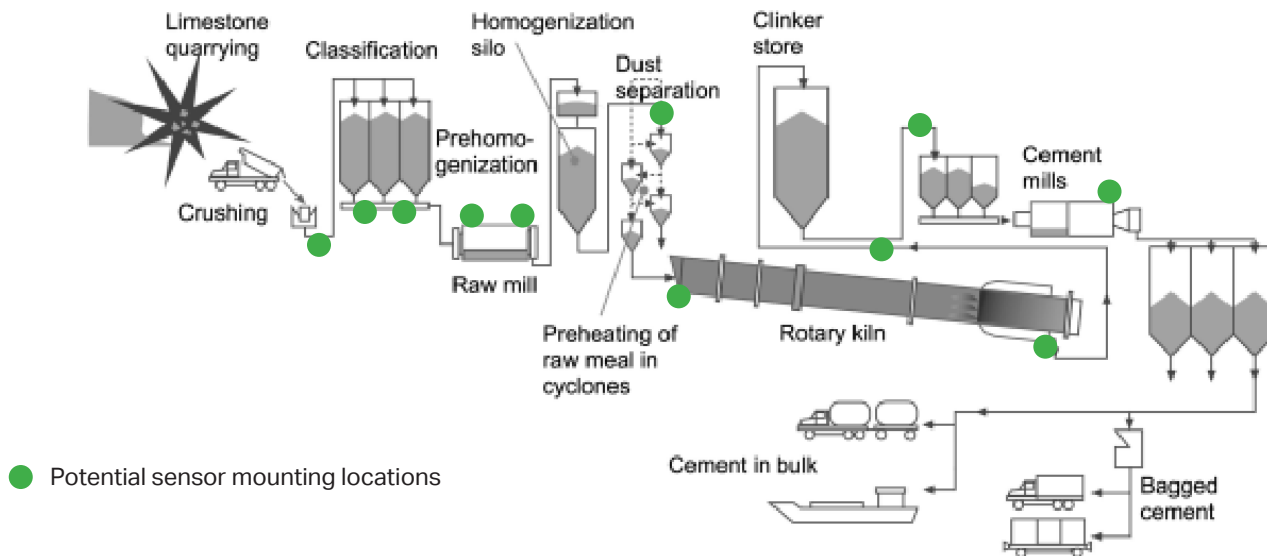
For customer service, call 1300-735-292
To email an order, ordersau@thermofisher.com
To order online: thermofisher.com

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

Data driven decision-making | Critical vibration measurement points | Thousands of sensors per plant



Key considerations, recommended solutions and products for cement factories

Consideration	Solution	Product
Imbalance on fans or motors	4-20 mA velocity sensors, standard accelerometers	PC420V, 786A
Contamination from dust and debris	Side exit sensors with strong hermetic seal, IP68 rated cable assemblies, stainless steel enclosures	787A, R6Q-J9T2AS, VLS junction boxes
High temperatures in kiln areas	High temperature accelerometers and cable/connector assemblies	HT780A, HT786A + R6Q-J9T2A
Belt drive misalignment	Tight tolerance accelerometers, intelligent vibration transmitter for dynamic analysis and connection to plant infrastructure	786A, 787A + iT300
Bearing damage or fatigue	Dynamic, high-frequency accelerometers and intelligent vibration transmitter for field-configurability and direct connection to control system	786A, 786A-I + iT300
Conveyors (running speed <3600 RPM)	4-20 mA sensors, high sensitivity accelerometers	PC420 sensors, 786-500, 787-500

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