

# Vibration monitoring for pulp & paper manufacturing

For over 60 years, Wilcoxon has provided solutions for the extreme industrial environments found on paper machines and has designed products that offer reliable long-term performance.

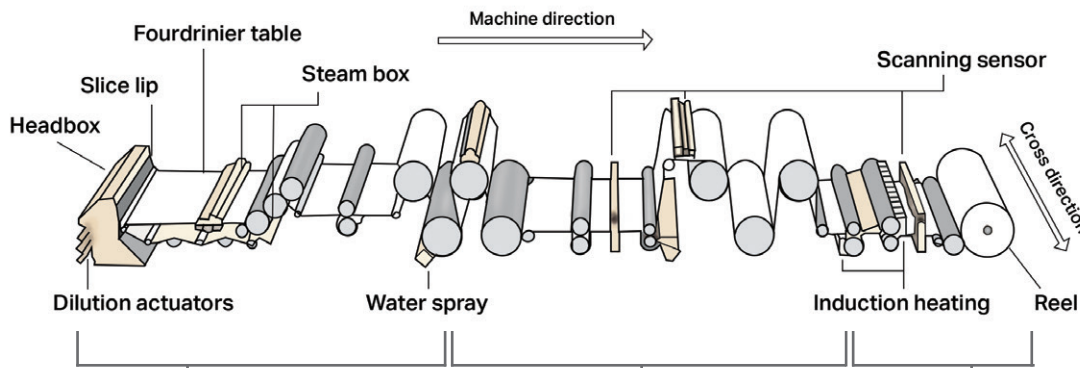


With the high number of rotating machines found in pulp and paper manufacturing facilities and their criticality to continuous operation, there's a huge potential benefit to employing vibration monitoring. High temperature accelerometers with reliable performance up to 165°C are ideal for use on rotating components in dryer sections, and rugged IP68 rated connectors and cables have the chemical and heat resistance needed throughout the plant.

## Considerations for the pulp & paper industry

- Consistently high operating temperatures cause sensor degradation
- High moisture levels damage electronics
- Caustic chemicals damage cabling
- Ensuring continuous operation, avoiding unplanned shutdowns

## Recommended products



Forming/press sections	Dryer section	Calendering section
Caustic chemicals, high moisture	High temperatures, sudden temperature spikes, high moisture	High moisture, high pressure
<ul style="list-style-type: none"> <li>• 786A/787A standard accelerometers</li> <li>• 786LF low frequency accelerometer</li> <li>• 6Q splash-proof connectors</li> <li>• J9T2A Teflon cables</li> </ul>	<ul style="list-style-type: none"> <li>• HT series high temperature accelerometers (FireFet)</li> <li>• 6Q splash-proof connectors</li> <li>• J9T2A Teflon cables</li> </ul>	<ul style="list-style-type: none"> <li>• 786A/787A accelerometers</li> <li>• 6Q/6W connectors</li> <li>• J9T2A/J9F Teflon cables</li> </ul>



## Authorised distributor

### In Australia:

For customer service, call 1300-735-292  
To email an order, [ordersau@thermofisher.com](mailto:ordersau@thermofisher.com)  
To order online: [thermofisher.com](http://thermofisher.com)

### In New Zealand:

For customer service, call 0800-933-966  
To email an order, [ordersnz@thermofisher.com](mailto:ordersnz@thermofisher.com)  
To order online: [thermofisher.com](http://thermofisher.com)