



Pharmaceutical applications

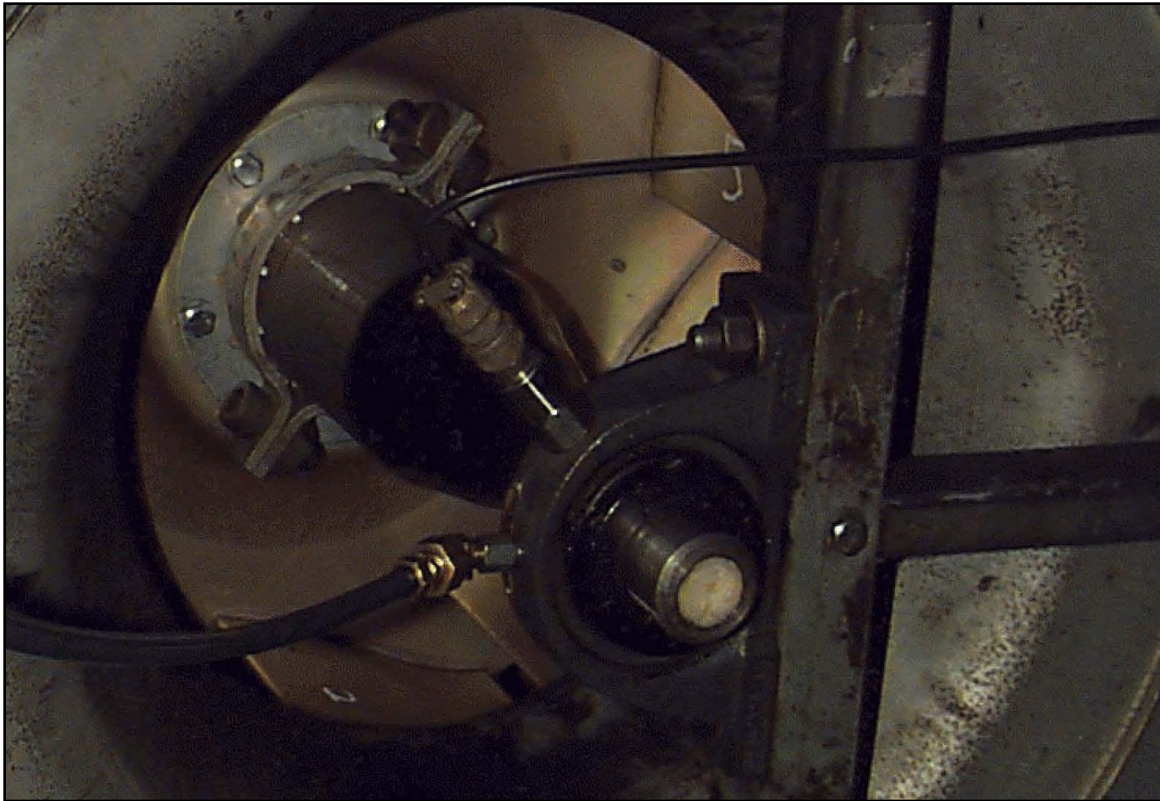
Wilcoxon
SENSING TECHNOLOGIES

Air handling

- Monitor
 - Flow turbulence (broadband energy, velocity)
 - Potential cavitation problems (broadband energy, acceleration)
 - Blade pass frequencies (BPF – discrete frequencies with harmonics)
 - Trim balance the rotors on the air handlers (1x running speed)



Air handler



Wilcoxon models used: 780A and 786A

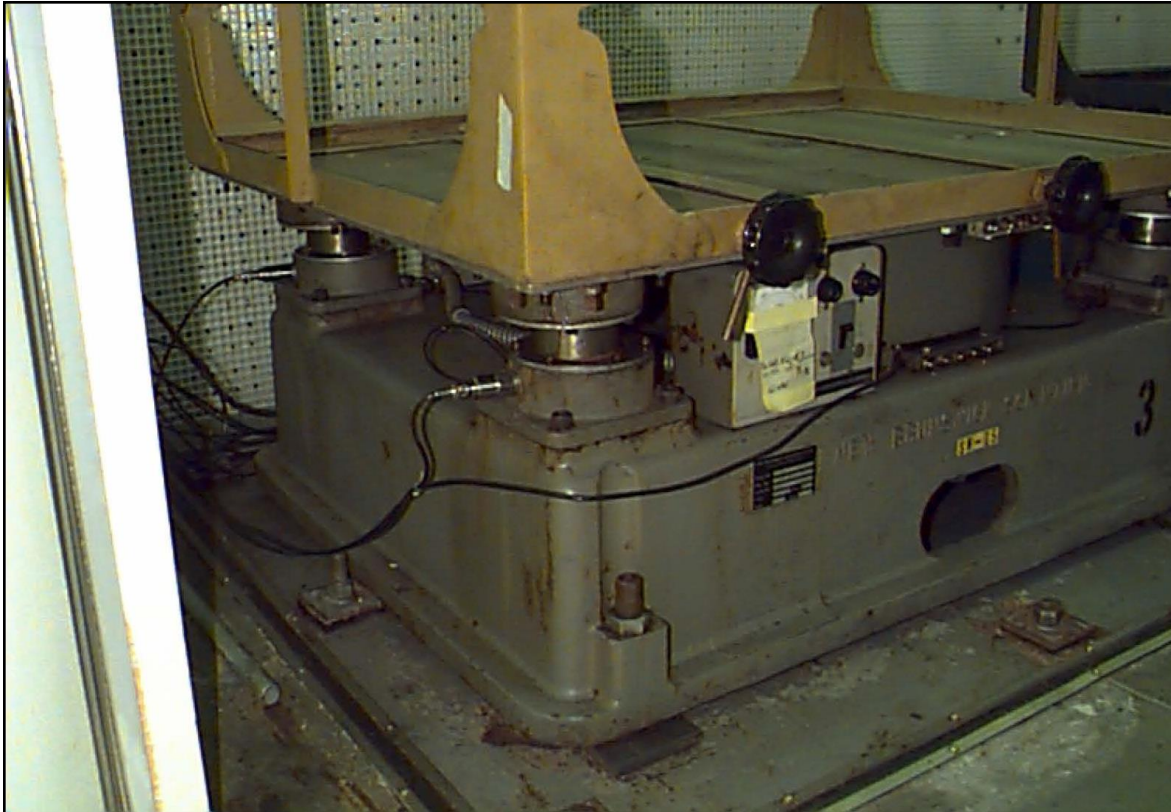


Chemical shaker

- Monitor
 - The bearings of a low vibration level chemical shaker (machinery maintenance – discrete frequencies)
 - The vibration level of the shaker (process measure – broadband energy)

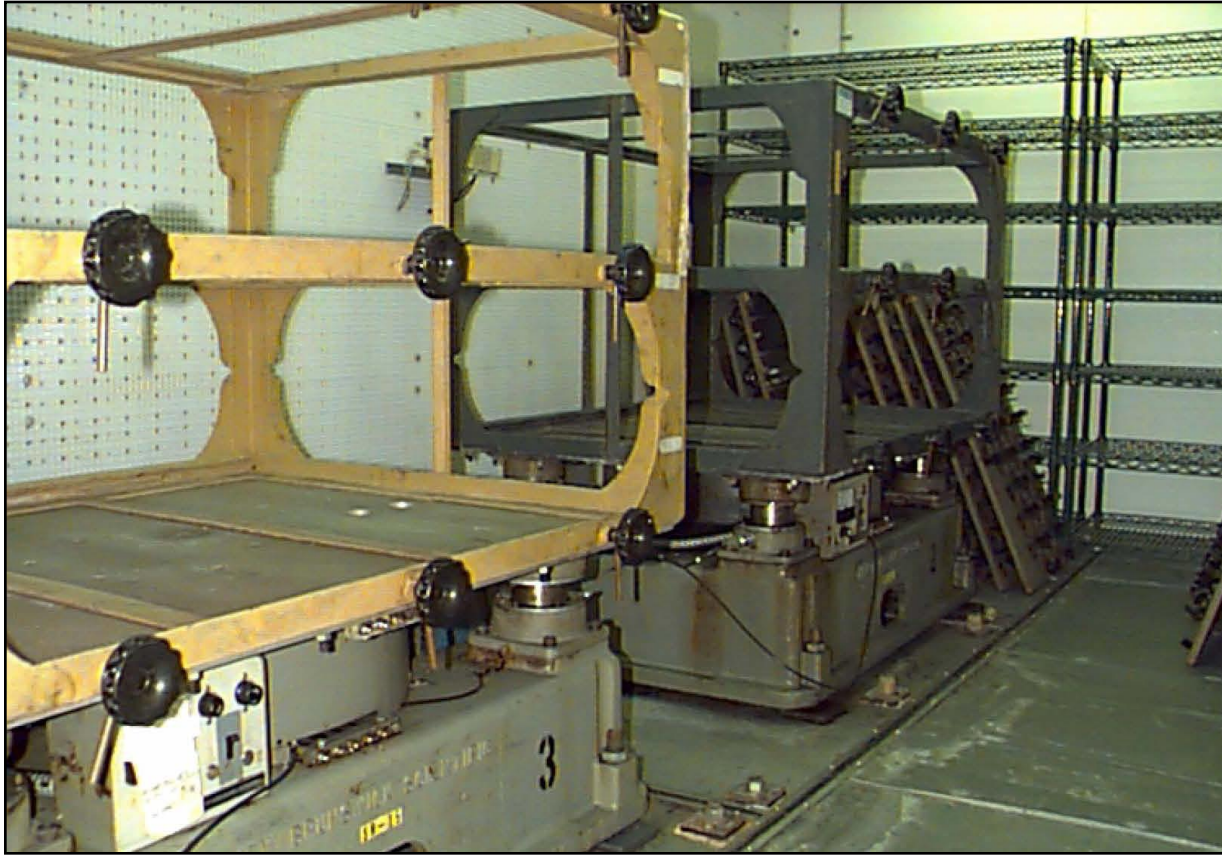


Chemical shaker



Wilcoxon model used: 786A

Chemical shaker



Chemical shaker

- Monitor
 - The bearings of a low vibration level chemical shaker (machinery maintenance – discrete frequencies)
 - The vibration level of the shaker (process measure – broadband energy)

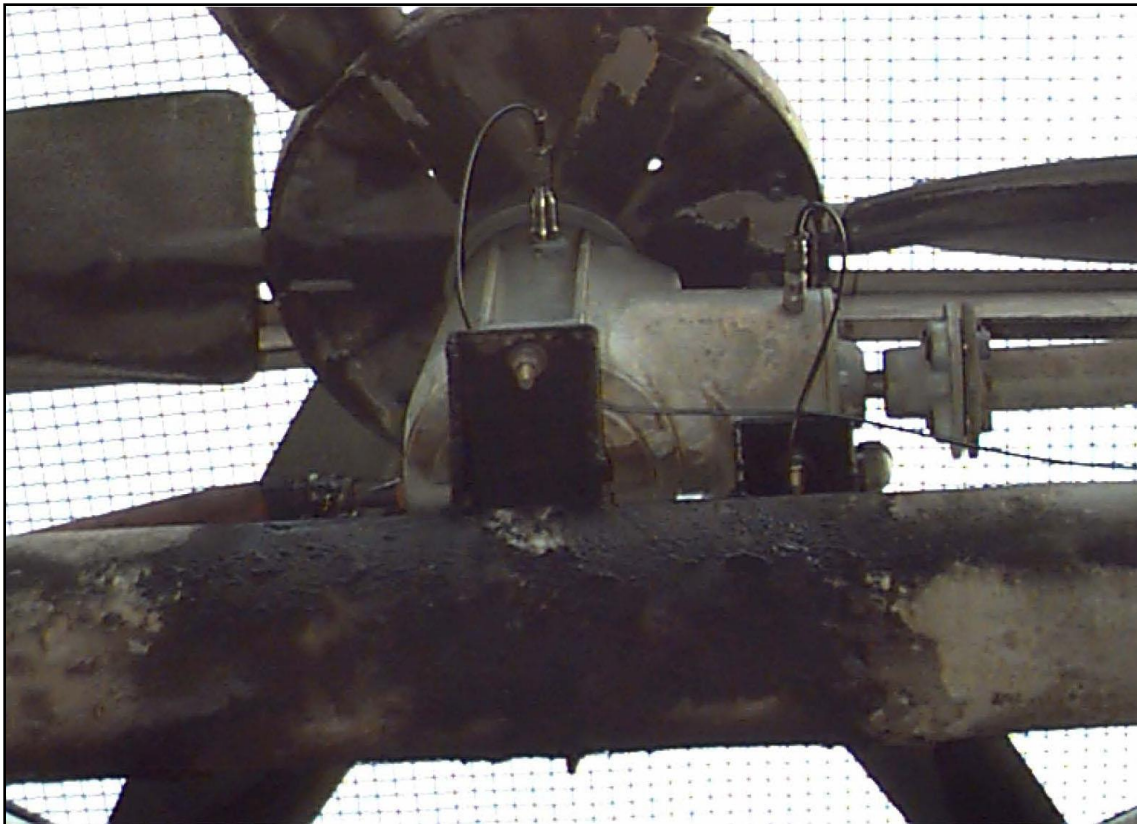


Cooling tower application

- Monitor
 - The shaft on a low speed fan for imbalance (gearbox output, <100 rpm)
 - Motors and gearboxes (1x, gearmesh and harmonics)
 - Blade imbalance (Low speed 1x)
 - Bearing wear (high frequency metal to metal contact)



Cooling tower fan



Wilcoxon model used: 793L and 786A

The information contained in this document is the property of Wilcoxon Sensing Technologies and is proprietary and/or copyright material. This information and this document may not be used or disclosed without the express authorization of Wilcoxon Sensing Technologies. Any unauthorized use or disclosure may be unlawful.

The information contained in this document may be subject to the provisions of the Export Administration Act of 1979 (50 USC 2401-2420), the Export Administration Regulations promulgated thereunder (15 CFR 730-774), and the International Traffic in Arms Regulations (22 CFR 120-130). The recipient acknowledges that these statutes and regulations impose restrictions on import, export, re-export and transfer to third countries of certain categories of data, technical services and information, and that licenses from the US Department of State and/or the US Department of Commerce may be required before such data, technical services and information can be disclosed. By accepting this document, the recipient agrees to comply with all applicable governmental regulations as they relate to the import, export and re-export of information.'

Authorised distributor

In Australia:

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

In New Zealand:

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

ThermoFisher
S C I E N T I F I C