



Machine tools and the automotive industry

Machine tool applications

- Transfer lines
- Grinding
- Milling
- Turning
- Broaching



Other automotive applications

- Casting
- Stamping
- Drawing (wire)
- Assembly



Machine tool – Spindle speed 10 to 300 Hz

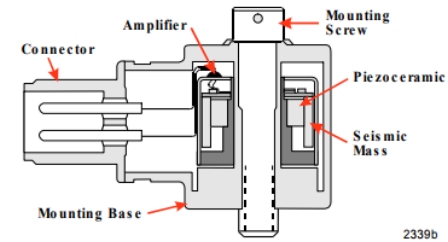
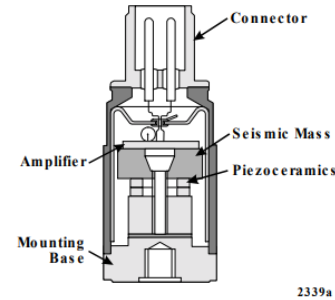
- Environment
 - Cutting fluid and cooling fluid spray, flying chips, electromagnetic interference
- Sensor requirements
 - Hermetic sealing
 - Splash proof / waterproof connectors
 - Teflon or Armored cabling
- Wilcoxon sensor recommendations
 - 793, 797, 786A – Industrial sensors
 - R6SLI, R6W – Sealed connectors
 - J9T2A cables



Slow speed machine tool – Spindle speed .5 to 10 Hz

- Sensor requirements
 - Very low noise
 - Very low strain sensitivity

- Wilcoxon sensor recommendations
 - 793L, 797L, 799LF – Industrial sensors
 - R6SLI, R6W – Sealed connectors
 - J9T2A cables



High speed machine tool – Spindle speed > 300 Hz


- Sensor requirements
 - Very high resonance
 - Low sensitivity
 - Overload protection

- Wilcoxon sensor recommendations
 - 732A
 - 732AT



Wilcoxon Sensing Technologies

For more information, please contact us:

 +1 (301) 330 8811

 info@wilcoxon.com

 www.wilcoxon.com

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