

Customer Focused, Quality Driven

Molecular Devices is committed to keeping our promises to customers and interested parties by:

- Continually improving our products, services, and processes
- Complying with applicable regulatory requirements
- Consistently meeting or exceeding our customer's expectations



Greg Milosevich
President, Molecular Devices

