



Customer Focused Quality Driven

Molecular Devices LLC 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 expectations



Mary Duseau

Mary Duseau | President | Molecular Devices

Contact Us

Phone: +1.800.635.5577
Web: www.moleculardevices.com
Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+91.73.8661.1198

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom





Global Environmental, Health and Safety Policy

Molecular Devices LLC is committed to conducting its business in an environmentally responsible manner, in compliance with all applicable environmental, health and safety (EHS) laws and regulations, and in a manner that promotes and protects the health and safety of our associates, contingent workers, visitors, contractors, customers, other business partners and members of our local communities worldwide.

We implement this policy by integrating environmental, health and safety requirements into operations through the Danaher Business System (DBS).

Molecular Devices operates in accordance with the following principles in conducting our business. These principles support our environmental, health and safety commitment:

- Compliance with all applicable environment, health and safety laws and regulations, including RoHS, REACH and applicable corporate, platform and operating company policies, standards and procedures.
- Associates are consulted with and encouraged to participate in the EHS management process, and feedback is solicited regarding ways to continually improve Molecular Devices' EHS program (through our annual associate engagement survey, Global EHS Community of Practice (which comprises EHS leaders from Danaher's operating companies around the world) and site-specific EHS Committees).
- Adoption of meaningful EHS priorities and objectives designed to drive continuous improvement, including with respect to prevention and reduction of workplace injuries and illnesses. Review of progress on a regular basis at multiple levels of the organization.
- Routine, formal evaluation of our compliance with applicable laws, regulations and policies, utilizing a combination of self-assessments and independent in-person audits.
- Integrity and accountability in personal conduct.
- Molecular Devices associates are expected to understand the environment, health and safety issues and risks associated with their jobs.
- Molecular Devices associates have a duty to act in an environmentally responsible and safe manner; this duty extends to fellow associates and contingent workers, visitors, contractors, customers, other business partners and members of our local communities.
- Commitment to continuous improvement by utilizing DBS-aligned EHS management systems and tools, including an EHS maturity assessment process and a suite of DBS EHS tools.
- Integration of sound environment, health and safety practices into applicable business functions to prioritize resources and risk mitigation efforts, with the goal of reducing incidents and preventing incidents that could result in life-threatening or life-altering outcomes.
- Consideration of environment, health and safety impacts in the course of developing new products or processes and in selecting production materials.
- Design, operation and maintenance of our facilities in a manner that minimizes greenhouse gas emissions and waste and protects water resources.
- Responsible use of materials, including, where feasible, the recycling and reuse of materials.
- EHS & sustainability due diligence assessments of all business acquisitions and before buying, leasing or selling property.
- Sensitivity to community concerns about environment, health, safety and sustainability issues.



Mary Duseau

Mary Duseau | President | Molecular Devices

Contact Us

Phone: +1.800.635.5577
 Web: www.moleculardevices.com
 Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+91.73.8661.1198

*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom

