

April 5th, 2010

RE: Molecular Devices Compliance with the Buy America Act

Dear Dr. Valued Customer:

Thank you for considering Molecular Devices, Inc. as your preferred partner for innovative life science research and drug discovery instruments, reagents and consumables.

Molecular Devices proudly maintains facilities in Sunnyvale, California and Downingtown, Pennsylvania. Our roots in the United States are deep. More than 20 years ago, we began offering our top-quality Microplate Readers to American researchers. Today, the **Molecular Devices division employs nearly 500 people in the United States**. Our employees work in over **15 states from Massachusetts to California**. Our sales representatives, field service engineers and field applications scientists live and work throughout the U.S. We have close partnerships with numerous U.S. firms from material/OEM component suppliers to ad agencies, raising the number of U.S. jobs that directly depend on Molecular Devices.

In addition, our **extensive portfolio of Life Science and Drug Discovery instruments, reagents and consumables comply with the “Buy America Act” for U.S. governmental purchases**. The Buy American Act (BAA - 41 U.S.C. § 10a–10d) requires the United States government to prefer U.S. or safe harbor country made products in its purchases.

All in all, Molecular Devices is proud of our contributions to, and participation in, the U.S. economy. We are deeply committed to helping the U.S. economy make a complete recovery. We celebrate the spirit and commitment to honesty, innovation and excellence that makes this country great. With over 100,000 products in operation in laboratories around the world, we look forward to working with you.

Regards,



Greg M. Milosevich
Vice President of Sales & Service, Americas
Molecular Devices
1311 Orleans Drive
Sunnyvale, CA 94089
www.moleculardevices.com
greg.milosevich@moldev.com
Mobile: 1-925-389-0771
Office: 1-408-548-6335