

1311 Orleans Drive
Sunnyvale, CA
94089-1136

Tel +1-800-635-5577
Fax +1-408-548-6461

Sales & Support

United States and Canada
1 800 635 5577

Brazil
+55 11 3616 6607

United Kingdom
+44 118 944 8000

Germany
+49 89/96 05 88 0

China
+86 10 6410 8669 (Beijing)
+86 21 6887 8820 Shanghai)

Japan
+81 6 6399 8211 (Osaka)
+81 3 5282 5261 (Tokyo)

South Korea
+82 2 3471 9531

Dear Molecular Devices Axon Instruments customer,

Thank you for your recent purchase. I am pleased to announce the availability of an important software update that addresses almost 20 issues reported by users like you. Please visit www.moleculardevices.com/pclamp10 to download pCLAMP™ or AxoScope software version 10.4.1.10.

I appreciate your loyalty to Axon Instruments and look forward to continuing to serve you in the future.

Best regards,

Jeffrey Tang, PhD
Product Manager of Axon Instruments