



May 28, 2019

To Whom it May Concern:

Starting year 2019, the manufacturing of the SpectraMax® M Series Multi-Mode Microplate Readers (M2/M2e, M3, M4, M5/M5e), FlexStation® 3 Multi-Mode Microplate Reader, Gemini™ XPS and EM, SpectraMax® L, Plus 384, 190, 340PC384, VersaMax™, and StakMax® Microplate Handling System will incorporate some cosmetic design updates to bring the product line into design alignment with our current brand guidelines.

These newly designed instruments will match the previous products in form, fit, and function. There are no changes to the internal workings of the instrument, this is merely an update to the colors of the LCD screen and/ or the colors of the outside chassis. The manufacturing process and verification testing remain identical and conform to our ISO 9001 standards.

We appreciate your loyalty to Molecular Devices readers and look forward to partnering with you now and in the future.

If you have any additional questions in this matter, please don't hesitate to reach out to contact our Technical Support team at support@moldev.com or 1-800-635-5577 in North America and techsupporteu@moldev.com or +44 118 944 8000 in Europe.

Sincerely,



Shreen Ibrahim
Director, Global Quality and EHS

5071967 Rev: B