

July 22, 2017

Statement Regarding the China RoHS Compliance of Molecular Devices Products

Please refer to Table 1 for the names and contents of the toxic or hazardous substances or elements contained in Molecular Devices products.

Table 1: Names and Contents of Toxic or Hazardous Substances or Elements

Part Name	Toxic or hazardous Substances and Elements					
	Lead (Pb)	Mercury (Hg)	Cadmium (Cd)	Hexavalent Chromium (Cr (VI))	PolyBrominated Biphenyls (PBB)	PolyBrominated Diphenyl Ethers (PBDE)
All Instruments	X	O	O	O	O	O

This table is prepared in accordance with the provisions of SJ/T 11364.

- O Indicates that said hazardous substance contained in all homogeneous materials for this part is below the limit requirement of GB/T 26572.
- X Indicates that said hazardous substance contained in at least one of the homogeneous materials used for this part is above the limit requirement of GB/T 26572.

Yours truly,




Raymond Mo
 Product Compliance Manager
 Molecular Devices, LLC

Legal Disclaimer:
 Molecular Devices LLC may make changes, improvements and/or modifications to the information provided at any time without notice.
 In NO EVENT shall Molecular Devices LLC, its employees, its officers, its affiliates, its suppliers or other third parties be liable for any damage, including without limitation special, indirect, or consequential damages (including, without limitation, damages for loss of profits, business interruption, or loss of information), arising out of the use, inability to use, or the results of use of information provided herein.