

Refer to the Bio-Plex hardware instruction manual for more details on the principles and procedures of calibration.

For technical service, call your local Bio-Rad office, or in the US, call 1-800-424-6563

Bio-Plex™ Calibration Kit Product Insert

For calibration of the Bio-Plex array reader

For research use only.
Not for diagnostic procedures.



Bio-Plex Array Reader Calibration

The Bio-Plex array reader requires daily calibration for optimal performance and reproducibility of results. This kit contains two calibrators (CAL1 and CAL2). These are microspheres with stable fluorescent intensities in the RP1, CL1 and CL2 wavelength ranges.

Target Values

Target values are printed on the label of each bottle. Enter these values into the calibration dialog in the Bio-Plex Manager™ software. See Bio-Plex Manager user guide for detailed instructions on entering target values and performing calibration.

Storage	2–8°C
Stability	6 months
Calibration frequency	Daily, or if the air temperature changes by more than $\pm 2^{\circ}\text{C}$ during the course of the day
CAL1	Calibrates the classification fluorescence detectors. Use the DD, CL1 and CL2 target values printed on the label
CAL2	Calibrates the reporter fluorescence detector. Use the RP1 target values printed on the label to calibrate the Bio-Plex system: Low RP1 target value: Bio-Plex cytokine assays High RP1: target value: Bio-Plex phosphoprotein assays