



Fluorescence Polarization SpectraMax Detection Cartridge

Fluorescence Polarization SpectraMax® Detection Cartridges offers great sensitivity for SpectraMax® i3 and Paradigm® Systems. The dual photomultiplier tube design on SpectraMax Paradigm accelerates the throughput 2X due to the direct excitation optical path.

