

# SpectraMax Paradigm



## Fluorescence Polarization Detection Cartridge



The **SpectraMax® Paradigm®** **Fluorescence Polarization Detection Cartridge** takes full advantage of the standard dual photomultiplier tube design of the SpectraMax Paradigm System to collect all the necessary data in a single read of a microplate. The resulting speed and performance are exceptional – whether reading in the top or bottom mode – due to the direct excitation optical path.

