

SpectraMax Paradigm



Tunable Wavelength Detection Cartridge



The **SpectraMax® Paradigm® TUNE Detection Cartridge** offers unmatched sensitivity in wavelength tunable detection. Designed to work with top or bottom read, it supports detection in:

- Fluorescence Intensity (FI)
- Time-Resolved Fluorescence (TRF)
- Luminescence (LUM)

