

Spark capabilities	GENOMICS AND PROTEOMICS	MICROBIOLOGY	CELL-BASED ASSAYS	DRUG DISCOVERY
Heating	●	●	●	●
Absorbance	●	●	●	●
Cuvette	●	●		●
Fluorescence filter top/bottom		●		
Fluorescence monochromator top/bottom	●		●	
Fluorescence Fusion Optics top/bottom				●
Fluorescence variable bandwidth			●	●
Fluorescence polarization				●
Fluorescence dichroic mirrors			●	●
Luminescence	●	●	●	●
Luminescence multicolor and scanning			●	●
Alpha Technology				●
Lid Lifter™		●	●	●
Cell counting and imaging		●	●	●
CO ₂ control				●
CO ₂ + O ₂ control		●	●	
Te-Cool™ cooling module				●
NanoQuant Plate™	●			