

Capabilities	SPARK CYTO 300	SPARK CYTO 400	SPARK CYTO 500	SPARK CYTO 600
Fluorescence Imaging	•	•	•	•
Brightfield imaging	•	•	•	•
Digital phase contrast imaging	•	•	•	•
Absorbance UV/vis monochromator - STD 384	•	•		
Absorbance UV/vis monochromator - ENH 1536			•	•
Fluorescence - STD 384	•			
Fluorescence - ENH 1536		•	•	•
Fluorescence filter top/bottom	•		•	
Fluorescence monochromator top/bottom		•		
Fluorescence Fusion Optics top/bottom				•
Fluorescence variable bandwidth		•		•
Fluorescence polarization		•	•	•
Fluorescence dichroic mirrors		•	•	•
Luminescence - STD 384 / multi-color and scanning	•	•		
Luminescence - ENH 1536 / multi-color and scanning			•	•
Alpha technology				•
Lid Lifter™	•	•	•	•
Heating	•	•	•	•
CO ₂ control	•	•	•	•
O ₂ control	•	•	•	•



Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170 **Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11

© 2020 Tecan trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com

