

Capabilities	SPARK CYTO 300	SPARK CYTO 400	SPARK CYTO 500	SPARK CYTO 600
Fluorescence imaging	•	•	•	•
Bright field imaging	•	•	•	•
Digital phase contrast imaging	•	•	•	•
Absorbance UV/vis monochromator	•	•	•	•
Fluorescence – standard	•			
Fluorescence – enhanced		•	•	•
Fluorescence filter top/bottom	•		•	•
Fluorescence monochromator top/bottom		•		•
Fluorescence variable bandwidth		•		•
Fluorescence polarization			•	•
Fluorescence dichroic mirrors		•	•	•
Luminescence	•	•	•	•
Luminescence multicolor and scanning			•	•
Alpha technology				•
Lid Lifter™	•	•	•	•
Heating	•	•	•	•
CO <sub>2</sub> and O <sub>2</sub> 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

© 2019 Tecan trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit [www.tecan.com](http://www.tecan.com)

