



Tecan: Spark/Spark Cyto Microplate Readers Configuration for Valita Aggregation Pure Assay

Parameter	Instrument Settings
Plate type	Corning 96-well Half Area Black Flat Bottom Non-binding, #3686
Mode	Fluorescence Polarization
Excitation	Filter (preferred) or monochromator
Excitation Wavelength	560nm
Excitation bandwidth	20nm
Emission	Filter (preferred) or monochromator
Emission Wavelength	600nm
Emission bandwidth	10nm
Flash Number	>= 100
Integration	40 μ s
Blank correction	None; 0mg/L standard should be labelled as standard and not blank
Gain	Optimized using most fluorescent well [0mg/L]
RFU (%)	70%
Z position	Optimized using most fluorescent well [0mg/L]
Mirror	Automatic
G-Factor	Default: 1.0 Manual adjustment may be required to set 0mg/L standard to approx. 70mP Increase G-factor to reduce mP, Decrease G-factor to increase mP

For research use only. Not for diagnostic purposes.

This configuration is for reference only and is not validated by Beckman Coulter. Beckman Coulter makes no warranties of any kind whatsoever expressed or implied, with respect to this configuration, including but not limited to warranties of fitness for a particular purpose or merchantability or that the configuration is non-infringing. All warranties are expressly disclaimed. Your use of the configuration is solely at your own risk, without recourse to Beckman Coulter.

© 2024 Beckman Coulter, Inc. All rights reserved. Valita, Valita Aggregation Pure and the ValitaCell logo are trademarks or registered trademarks of ValitaCell Ltd. in the United States and other countries. ValitaCell is a Beckman Coulter Company. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

2024-GBL-EN-104996-v1