



Agilent: Biotek Synergy Neo 2 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 Polarisation
PMT	Synergy Neo 2S: Single PMT Synergy Neo 2: Dual PMT
Filter Cube	Synergy Neo 2S: Single PMT Red Cube (#120: EX 530/25, EM 590/35, DM570) Synergy Neo 2: Dual FP Red Cube (#62: EX 530/25, EM 590/35, DM570) Additional upper cube (#4) also required
Read Speed	Normal
Delay after plate movement	0 msec
Measurement per data point	≥ 100
Gain	'Options' > 'Automatic Gain Adjustment' > 'Scale to High wells' and 10% target value. Adjustment using most fluorescent well [0mg/L]
Read Height	Options > automatic adjustment using most fluorescent well [0mg/L]
Dynamic Range	'Standard' (0 to 99,999)
Lamp Energy	High
G-Factor	Default 1 can be used, but manual adjustment preferred
Blank correction	None; 0mg/L standard should be labelled as standard and not blank

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-104994-v1