

Product Number:	101351
Lot Number:	504115
Expiration:	28-Jun-2024

Component	Part Number	Lot Number	Control Number	Concentration
488-dyed Singleplex Beads ²	104007	329215	N/A	1.23 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	319815	319815A	
Bead Diluent	100458	333912	N/A	
SBG Dilution Buffer	100376	330321	N/A	
SBG Concentrate	103397	319210	319210A	50 nM
Bead Conjugation Buffer	100453	402307	N/A	
Bead Wash Buffer	101352	318108	318108A	
Bead Blocking Buffer	100457	402510	N/A	
Biotinylation Reaction Buffer	100460	334009	N/A	
EDC, 10mg	100022	RM2023060000068	N/A	
EZ-Link NHS-PEG4 Biotin	100348	RM2024010000049	N/A	
RGP ¹	103159	N/A ¹	N/A ¹	
PSA Capture Antibody	100463	332606	N/A	5.1 mg/mL
PSA Detector Antibody	100464	402605	N/A	1.1 mg/mL
Calibrator A-H	100758 - 100765	305311	N/A	
Control 1	100768	305312	N/A	
Control 2	100769	305312	N/A	

¹ Reagents are not Kit Lot Specific

² Assay Beads provided are conjugated with 488 nm fluorescent dye.

Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Spec. Range (pg/mL)	
Control 1	305312	2.048 - 3.072	
Control 2	305312	41.137 - 65.144	

Calibrator Levels (pg/mL)	
A	0.000
B	0.077
C	0.246
D	0.873
E	2.61
F	8.43
G	25.8
H	64.4

Review/Approval

ROBERT SOMOGTI

QC MANAGER

[Signature]

05 FEB 2024

Name

Title

Signature/Date