

Product Number:	101351
Lot Number:	504127
Expiration:	28-Jun-2024

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads <sup>2</sup>	104007	329215	1.23 x 10 <sup>9</sup> beads/mL
Detector / Sample Diluent	101353	322620	
Bead Diluent	100458	333912	
SBG Dilution Buffer	100376	330321	
SBG Concentrate	103397	330603	50 nM
Bead Conjugation Buffer	100453	328909	
Bead Wash Buffer	101352	323415	
Bead Blocking Buffer	100457	320109	
Biotinylation Reaction Buffer	100460	334009	
EDC, 10mg	100022	RM2023060000068	
EZ-Link NHS-PEG4 Biotin	100348	RM2024010000049	
RGP <sup>1</sup>	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	332606	5.1 mg/mL
PSA Detector Antibody	100464	330605	1.1 mg/mL
Calibrator A-H	100758 - 100765	305311	
Control 1	100768	305312	
Control 2	100769	305312	

<sup>1</sup> Reagents are not Kit Lot Specific


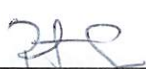
<sup>2</sup> Assay Beads provided are conjugated with 488 nm fluorescent dye.

Release Materials*		
	Lot	Spec. Range (pg/mL)
Control 1	305312	2.048 - 3.072
Control 2	305312	41.137 - 65.144

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

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



  
 Name: ROBERT SOMOGYI Title: QC MANAGER Signature/Date: 16 FEB 2024