

<b>Product Number:</b>	101351
<b>Lot Number:</b>	504341
<b>Expiration:</b>	22-Feb-2025

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads <sup>2</sup>	104007	403913	1.14 x 10 <sup>9</sup> beads/mL
Detector / Sample Diluent	101353	404108	
Bead Diluent	100458	412312	
SBG Dilution Buffer	100376	417210	
SBG Concentrate	103397	420404	50 nM
Bead Conjugation Buffer	100453	416608	
Bead Wash Buffer	101352	415103	
Bead Blocking Buffer	100457	411728	
Biotinylation Reaction Buffer	100460	405509	
EDC, 10mg	100022	RM2024070000011	
EZ-Link NHS-PEG4 Biotin	100348	RM2024060000021	
RGP <sup>1</sup>	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	414114	5.1 mg/mL
PSA Detector Antibody	100464	406102	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*			* 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	2.05 - 3.07	
Control 2	305312	41.1 - 65.1	

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 SONOGYI QC MANAGER [Signature] 05 AUG 2024

Name Title Signature/Date