

Product Number:	101351
Lot Number:	504154
Expiration:	30-Jan-2025

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	334113	1.18 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	402218	
Bead Diluent	100458	402924	
SBG Dilution Buffer	100376	402514	
SBG Concentrate	103397	401702	50 nM
Bead Conjugation Buffer	100453	402307	
Bead Wash Buffer	101352	403115	
Bead Blocking Buffer	100457	402510	
Biotinylation Reaction Buffer	100460	401708	
EDC, 10mg	100022	RM2024020000041	
EZ-Link NHS-PEG4 Biotin	100348	RM2024010000049	
RGP ¹	103159	N/A ¹	
PSA Capture Antibody	100463	402922	5.1 mg/mL
PSA Detector Antibody	100464	402605	1.1 mg/mL
Calibrator A-H	100758 - 100765	305311	
Control 1	100768	305312	
Control 2	100769	305312	

¹ Reagents are not Kit Lot Specific



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

Release Materials*		
	Lot	Spec. Range (pg/mL)
Control 1	100768	2.048 - 3.072
Control 2	100769	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: [Signature] 16 FEB 2024