

Product Number:	101351
Lot Number:	504272
Expiration:	22-Feb-2025

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	403913	1.14 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	412313	
Bead Diluent	100458	412312	
SBG Dilution Buffer	100376	407412	
SBG Concentrate	103397	408804	50 nM
Bead Conjugation Buffer	100453	409213	
Bead Wash Buffer	101352	403115	
Bead Blocking Buffer	100457	411728	
Biotinylation Reaction Buffer	100460	405509	
EDC, 10mg	100022	RM2024050000045	
EZ-Link NHS-PEG4 Biotin	100348	RM2023030000060	
RGP ¹	103159	N/A ¹	
PSA Capture Antibody	100463	402922	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	

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	305312	2.05 - 3.07
Control 2	305312	41.1 - 65.1

* 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

Robert Somogyi

QC Manager

 24 MAY 2024

Name

Title

Signature/Date