

Product Number:	101351
Lot Number:	503160
Expiration:	15-Apr-2022

Component	Part Number	Lot Number	Concentration
Beads	103611	115514	3.18 X 10 ⁹ beads/mL
Detector/Sample Diluent	101353	122113	
Bead Diluent	100458	118710	
SBG Dilution Buffer	100376	123912	
SBG Concentrate	103397	120031	50nM
Bead Conjugation Buffer	100453	120810	
Bead Wash Buffer	101352	117225	
Bead Blocking Buffer	100457	110524	
Biotinylation Reaction Buffer	100460	117226	
EDC, 10mg	100022	RM2021050000061	
EZ-Link NHS-PEG4 Biotin	100348	RM2021070000039	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	122803	5.1 mg/ml
PSA Detector Antibody	100464	122804	1.0 mg/ml
Calibrator A-H	100758-100765	121603	
Control 1	100768	121604	
Control 2	100769	121604	

¹ RGP is not Kit Lot specific.

	Mean Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	1.69-3.70	
Control 2	42.7-64.0	

Calibrator Levels (pg/mL)

- A 0.000
- B 0.043
- C 0.237
- D 0.937
- E 2.62
- F 8.76
- G 26.2
- H 61.6

Review/Approval

Jungil Park QC Supervisor [Signature] 28 Sep 2021
Name Title Signature/Date