

Product Number:	101351
Lot Number:	502850
Expiration:	28-Aug-2021

Component	Part Number	Lot Number	Concentration
Beads	103611	102533	2.98X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	105014	
Bead Diluent	100458	105417	
SBG Dilution Buffer	100376	108322	
SBG Concentrate	103397	104730	50nM
Bead Conjugation Buffer	100453	100708	
Bead Wash Buffer	101352	031122	
Bead Blocking Buffer	100457	024110	
Biotinylation Reaction Buffer	100460	100822	
EDC, 10mg	100022	RM2020120000032	
EZ-Link NHS-PEG4 Biotin	100348	RM2020120000031	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	105804	5.2 mg/ml
PSA Detector Antibody	100464	108106	1.0 mg/mL
Calibrator A-H	100758-100765	027208	
Control 1	100768	027209	
Control 2	100769	027209	

<sup>1</sup> RGP is not Kit Lot specific.

	Mean Range (pg/mL)	
Control 1	2.25-3.38	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 2	47.1-70.6	

**Calibrator Levels (pg/mL)**

A	0.000
B	0.095
C	0.281
D	0.943
E	2.95
F	9.81
G	28.5
H	73.6

Review/Approval

*Shivani Goel*  
Name

QC Manager  
Title

*Shivani Goel 07Apr 2021*  
Signature/Date