

Product Number:	101351
Lot Number:	502854
Expiration:	28-Aug-2021

Component	Part Number	Lot Number	Concentration
Beads	103611	102534	3.05X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	105014	
Bead Diluent	100458	105414	
SBG Dilution Buffer	100376	104729	
SBG Concentrate	103397	101913	50nM
Bead Conjugation Buffer	100453	100706	
Bead Wash Buffer	101352	031132	
Bead Blocking Buffer	100457	024112	
Biotinylation Reaction Buffer	100460	100821	
EDC, 10mg	100022	RM2020120000032	
EZ-Link NHS-PEG4 Biotin	100348	RM2020120000031	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	103512	5.1 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

*Jumpil Park*  
Name

QC Supervisor  
Title

*[Signature]* 08 Apr 2021  
Signature/Date