

Product Number:	101351
Lot Number:	503098
Expiration:	29-Mar-2022

Component	Part Number	Lot Number	Concentration
Beads	103611	115512	2.61X10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	119417	
Bead Diluent	100458	118710	
SBG Dilution Buffer	100376	121714	
SBG Concentrate	103397	119416	50nM
Bead Conjugation Buffer	100453	108831	
Bead Wash Buffer	101352	117225	
Bead Blocking Buffer	100457	110522	
Biotinylation Reaction Buffer	100460	117224	
EDC, 10mg	100022	RM2021050000061	
EZ-Link NHS-PEG4 Biotin	100348	RM2021070000039	
RGP	103159	N/A <sup>1</sup>	
PSA Capture Antibody	100463	119303	5.1 mg/ml
PSA Detector Antibody	100464	119304	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

*Wankai Tsai*

Name

QC

Title

*[Signature]* 11 Aug 2021

Signature/Date