

Product Number:	101351
Lot Number:	502505
Expiration:	15-Jan-2021

Component	Part Number	Lot Number	Concentration
<b>Beads</b>	103611	005515	3.07X10 <sup>9</sup> beads/mL
<b>Detector/Sample Diluent</b>	101353	009507	
<b>Bead Diluent</b>	100458	001517	
<b>SBG Dilution Buffer</b>	100376	005223	
<b>SBG Concentrate</b>	103397	002715	50nM
<b>Bead Conjugation Buffer</b>	100453	011707	
<b>Bead Wash Buffer</b>	101352	014004	
<b>Bead Blocking Buffer</b>	100457	002713	
<b>Biotinylation Reaction Buffer</b>	100460	001504	
<b>EDC, 10mg</b>	100022	RM2020020000021	
<b>EZ-Link NHS-PEG4 Biotin</b>	100348	RM2020040000009	
<b>RGP</b>	103159	N/A <sup>1</sup>	
<b>PSA Capture Antibody</b>	100463	006901	5.0 mg/ml
<b>PSA Detector Antibody</b>	100464	935009	1.1 mg/mL
<b>Calibrator A-H</b>	100758-100765	921201	
<b>Control 1</b>	100768	921202	
<b>Control 2</b>	100769	921202	

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

	Mean Range (pg/mL)	
Control 1	2.16-3.24	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 2	51.2-76.8	

**Calibrator Levels (pg/mL)**

A	0.000
B	0.084
C	0.252
D	0.899
E	2.82
F	9.94
G	30.0
H	76.3

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel 20 Aug 2020*

Name

Title

Signature/Date