

Product Number:	101351
Lot Number:	502586
Expiration:	18-Mar-2021

Component	Part Number	Lot Number	Concentration
Beads	103611	024812	3.17X10 ⁹ beads/mL
Detector/Sample Diluent	101353	028406	
Bead Diluent	100458	022412	
SBG Dilution Buffer	100376	026619	
SBG Concentrate	103397	019712	50nM
Bead Conjugation Buffer	100453	028402	
Bead Wash Buffer	101352	025907	
Bead Blocking Buffer	100457	024107	
Biotinylation Reaction Buffer	100460	025910	
EDC, 10mg	100022	RM2020030000046	
EZ-Link NHS-PEG4 Biotin	100348	RM2020040000009	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	019803	5.2 mg/ml
PSA Detector Antibody	100464	028405	1.1 mg/mL
Calibrator A-H	100758-100765	027208	
Control 1	100768	027209	
Control 2	100769	027209	

¹ 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	QC Manager	<i>Shivani Goel 23oct 2020</i>
	Name	Title	Signature/Date