

Product Number:	101351
Lot Number:	503054
Expiration:	24-Mar-2022

Component	Part Number	Lot Number	Concentration
Beads	103611	108830	2.60X10 ⁹ beads/mL
Detector/Sample Diluent	101353	111312	
Bead Diluent	100458	110117	
SBG Dilution Buffer	100376	115221	
SBG Concentrate	103397	108324	50nM
Bead Conjugation Buffer	100453	108831	
Bead Wash Buffer	101352	111608	
Bead Blocking Buffer	100457	109014	
Biotinylation Reaction Buffer	100460	117224	
EDC, 10mg	100022	RM2021050000020	
EZ-Link NHS-PEG4 Biotin	100348	RM2021070000039	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	105804	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	

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
Name

QC
Title

Shivani Goel 20/01/2021
Signature/Date