

Product Number:	101351
Lot Number:	503208
Expiration:	15-Apr-2022

Component	Part Number	Lot Number	Concentration
Beads	103611	108830	2.60 X 10 ⁹ beads/mL
Detector/Sample Diluent	101353	125711	50nM
Bead Diluent	100458	122710	
SBG Dilution Buffer	100376	128906	
SBG Concentrate	103397	120035	
Bead Conjugation Buffer	100453	120810	
Bead Wash Buffer	101352	122509	
Bead Blocking Buffer	100457	110524	
Biotinylation Reaction Buffer	100460	117226	
EDC, 10mg	100022	RM2021090000073	
EZ-Link NHS-PEG4 Biotin	100348	RM2021070000039	
RGP	103159	N/A ¹	5.1 mg/ml
PSA Capture Antibody	100463	119303	
PSA Detector Antibody	100464	108106	1.0 mg/ml
Calibrator A-H	100758-100765	121603	
Control 1	100768	121604	
Control 2	100769	121604	

¹ RGP is not Kit Lot specific.

	Mean Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	1.69-3.70	
Control 2	42.7-64.0	

Calibrator Levels (pg/mL)

A	0.000
B	0.043
C	0.237
D	0.937
E	2.62
F	8.76
G	26.2
H	61.6

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 04Nov2021
Signature/Date