

Product Number:	101351
Lot Number:	502269
Expiration:	16-Oct-2020

Component	Part Number	Lot Number	Concentration
Beads	103206	001510	3.24X10 ⁹ beads/mL
Detector/Sample Diluent	101353	001317	
Bead Diluent	100458	001517	
SBG Dilution Buffer	100376	001516	
SBG Concentrate	103397	935218	50nM
Bead Conjugation Buffer	100453	932512	
Bead Wash Buffer	101352	928911	
Bead Blocking Buffer	100457	002713	
Biotinylation Reaction Buffer	100460	935217	
EDC, 10mg	100022	RM2019120000025	
EZ-Link NHS-PEG4 Biotin	100348	RM2019120000049	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	903614	5.0 mg/ml
PSA Detector Antibody	100464	935009	1.1 mg/mL
Calibrator A-H	100758-100765	921201	
Controls 1 and 2	100768-100769	921202	

¹ 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

Patrizia Stadler Assoc. Director, QC *P. Stadler* 16 MAR 2020
 Name Title Signature/Date