

Product Number:	101351
Lot Number:	502043
Expiration:	5-Feb-2020

Component	Part Number	Lot Number	Concentration
Beads	103206	823620	3.41X10 ⁹ beads/mL
Detector/Sample Diluent	101353	924008	
Bead Diluent	100458	917022	
SBG Dilution Buffer	100376	921117	
SBG Concentrate	103397	921713	50nM
Bead Conjugation Buffer	100453	919112	
Bead Wash Buffer	101352	903602	
Bead Blocking Buffer	100457	920006	
Biotinylation Reaction Buffer	100460	917619	
EDC, 10mg	100022	RM2019070000066	
EZ-Link NHS-PEG4 Biotin	100348	RM2019090000016	
RGP	103159	N/A ¹	
PSA Capture Antibody	100463	925412	5.0 mg/mL
PSA Detector Antibody	100464	919203	1.0 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 27 Sep 2019*

Name	Title	Signature/Date
------	-------	----------------