

Product Number:	101351
Lot Number:	504010
Expiration:	24-Jan-2024

Component	Part Number	Lot Number	Control Number	Concentration
488-dyed Singleplex Beads ²	104007	307212	N/A	1.32 x 10 ⁹ beads/mL
Detector/Sample Diluent	101353	303216	N/A	
Bead Diluent	100458	312905	312905A	
SBG Dilution Buffer	100376	327510	N/A	
SBG Concentrate	103397	315713	315713A	50nM
Bead Conjugation Buffer	100453	306010	N/A	
Bead Wash Buffer	101352	302410	N/A	
Bead Blocking Buffer	100457	304807	N/A	
Biotinylation Reaction Buffer	100460	320502	N/A	
EDC, 10mg	100022	RM2023050000024	N/A	
EZ-Link NHS-PEG4 Biotin	100348	RM2023020000034	N/A	
RGP ¹	103159	N/A	N/A	
PSA Capture Antibody	100463	325002	N/A	5.1 mg/ml
PSA Detector Antibody	100464	320612	N/A	1.1 mg/ml
Calibrator A-H	100758-100765	305311	N/A	
Control 1	100768	305312	N/A	
Control 2	100769	305312	N/A	

¹RGP is not Kit Lot specific.

²Assay Beads provided are conjugated with 488 nm fluorescent dye.

	Spec. Range (pg/mL)	
Control 1	2.048 - 3.072	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 2	41.137 - 65.144	

Calibrator Levels (pg/mL)

- A 0.00
- B 0.077
- C 0.246
- D 0.873
- E 2.61
- F 8.43
- G 25.8
- H 64.4

Review/Approval: Sonia Khillan Name QC manager Title Sonia Khillan Signature/Date 16 OCT 2023