

Product Number:	101351
Lot Number:	503975
Expiration:	26-Oct-2023

Component	Part Number	Lot Number	Control Number	Concentration
488-dyed Singleplex Beads <sup>2</sup>	104007	312408	312408A	1.27 x 10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	319815	319815A	
Bead Diluent	100458	318009	318009A	
SBG Dilution Buffer	100376	316508	316508A	
SBG Concentrate	103397	315712	315712A	50nM
Bead Conjugation Buffer	100453	319116	319116A	
Bead Wash Buffer	101352	318108	318108A	
Bead Blocking Buffer	100457	320106	N/A	
Biotinylation Reaction Buffer	100460	320502	N/A	
EDC, 10mg	100022	RM2023050000024	N/A	
EZ-Link NHS-PEG4 Biotin	100348	RM2023040000067	N/A	
RGP <sup>1</sup>	103159	N/A	N/A	
PSA Capture Antibody	100463	303212	N/A	5.1 mg/ml
PSA Detector Antibody	100464	303213	N/A	1.1 mg/ml
Calibrator A-H	100758-100765	213502	N/A	
Control 1	100768	213503	N/A	
Control 2	100769	213503	N/A	

<sup>1</sup>RGP is not Kit Lot specific.

<sup>2</sup>Assay Beads provided are conjugated with 488 nm fluorescent dye.

For more information, please visit

<https://go.quanterix.com/HomebrewBeadUpdate>



	Spec. Range (pg/mL)	* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
Control 1	2.09 - 3.14	
Control 2	45.9 - 68.9	

Calibrator Levels (pg/mL)	
A	0.000
B	0.090
C	0.260
D	0.878
E	2.71
F	9.56
G	26.9
H	70.0

Review/Approval: Sonia Khillan Name: Sonia Khillan Title: QC Manager Signature/Date: Sonia Khillan 11/SEP/2023