

Product Number:	101351
Lot Number:	503983
Expiration:	9-Mar-2024

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	322620	N/A	
Bead Diluent	100458	306809	N/A	
SBG Dilution Buffer	100376	308609	N/A	
SBG Concentrate	103397	315713	315713A	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	RM2023060000067	N/A	
EZ-Link NHS-PEG4 Biotin	100348	RM2023040000067	N/A	
RGP <sup>1</sup>	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	

<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	2.048 - 3.072	
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: Sonia Khillan Title: QC Manager Signature/Date: Sonia Khillan 20SEP2023