

Product Number:	101351
Lot Number:	503993
Expiration:	1-Feb-2024

Component	Part Number	Lot Number	Control Number	Concentration
488-dyed Singleplex Beads ²	104007	312408	312408A	1.27 x 10 ⁹ beads/mL
Detector/Sample Diluent	101353	313512	313512A	
Bead Diluent	100458	318009	318009A	
SBG Dilution Buffer	100376	308609	N/A	
SBG Concentrate	103397	315713	315713A	50nM
Bead Conjugation Buffer	100453	303210	N/A	
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 ¹	103159	N/A	N/A	
PSA Capture Antibody	100463	325002	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	

¹RGP is not Kit Lot specific.

²Assay 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, QC Manager Title, Sonia Khillan 02OCT2023 Signature/Date