

Product Number:	101351
Lot Number:	503751
Expiration:	13-Apr-2023

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads <sup>2</sup>	104007	230417	1.25 x 10 <sup>9</sup> beads/mL
Detector/Sample Diluent	101353	231914	
Bead Diluent	100458	232109	
SBG Dilution Buffer	100376	231814	
SBG Concentrate	103397	222904	50nM
Bead Conjugation Buffer	100453	217503	
Bead Wash Buffer	101352	210923	
Bead Blocking Buffer	100457	210318	
Biotinylation Reaction Buffer	100460	215211	
EDC, 10mg	100022	RM2022100000009	
EZ-Link NHS-PEG4 Biotin	100348	XH359829	
RGP <sup>1</sup>	103159	N/A	
PSA Capture Antibody	100463	204106	5.1 mg/ml
PSA Detector Antibody	100464	225612	1.0 mg/ml
Calibrator A-H	100758-100765	121603	
Control 1	100768	121604	
Control 2	100769	121604	

<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	1.70 - 3.55	
Control 2	43.1 - 64.6	

Calibrator Levels (pg/mL)
A 0
B 0.043
C 0.237
D 0.937
E 2.62
F 8.76
G 26.2
H 61.6

Review/Approval

Sonia Khillan  
Name

QC  
Title

Sonia Khillan 05DEC2022  
Signature/Date