

Product Number:	101351
Lot Number:	503995
Expiration:	17-Feb-2024

Component	Part Number	Lot Number	Control Number	Concentration
488-dyed Singleplex Beads ²	104007	220210	N/A	1.34 x 10 ⁹ beads/mL
Detector/Sample Diluent	101353	308307	N/A	
Bead Diluent	100458	318009	318009A	
SBG Dilution Buffer	100376	319209	319209A	
SBG Concentrate	103397	315713	315713A	50nM
Bead Conjugation Buffer	100453	319116	319116A	
Bead Wash Buffer	101352	318108	318108A	
Bead Blocking Buffer	100457	304807	N/A	
Biotinylation Reaction Buffer	100460	306812	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	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.

For more information, please visit

<https://go.quoterix.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

QC Manager

Title

Sonia Khillan 04/06/2023

Signature/Date