

Product Number:	101351
Lot Number:	504626
Expiration:	30-Apr-2026

Component	Part Number	Lot Number	Control Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	513911	N/A	1.20 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	511806	N/A	50 nM
Bead Diluent	100458	514005	N/A	
SBG Dilution Buffer	100376	510511	510511A	
SBG Concentrate	103397	506201	506201A	
Bead Conjugation Buffer	100453	509411	N/A	
Bead Wash Buffer	101352	509410	N/A	
Bead Blocking Buffer	100457	513908	N/A	
Biotinylation Reaction Buffer	100460	510809	N/A	
EDC, 10mg	100022	RM2025040000039	N/A	
EZ-Link NHS-PEG4 Biotin	100348	RM2025050000023	N/A	
RGP ¹	103159	N/A	N/A ¹	

¹ Reagents are not Kit Lot Specific

² Assay Beads provided are conjugated with 488 nm fluorescent dye.

Review/Approval

Dhruv Patel

DOC Control/II, QA

02 Jun 2025

Name

Title

Signature/Date

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, 01821

techsupport@quanterix.com

©2025 Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures.

Release

COA-0057 02