

Product Number:	101351
Lot Number:	504700
Expiration:	30-Apr-2026

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	519910	1.26 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	511808	50 nM
Bead Diluent	100458	512007	
SBG Dilution Buffer	100376	515316	
SBG Concentrate	103397	515406	
Bead Conjugation Buffer	100453	513606	
Bead Wash Buffer	101352	521010	
Bead Blocking Buffer	100457	521011	
Biotinylation Reaction Buffer	100460	510810	
EDC, 10mg	100022	RM2025040000039	
EZ-Link NHS-PEG4 Biotin	100348	RM2025050000023	
RGP ¹	103159	N/A ¹	

¹ Reagents are not Kit Lot Specific

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

Review/Approval

Dharmi Patel QA Specialist II, QA

DR / 03 Sep 2025.

Name

Title

Signature/Date