

Product Number:	101351
Lot Number:	504739
Expiration:	16-May-2026

Component	Part Number	Lot Number	Concentration
488-dyed Singleplex Beads ²	104007	520513	1.29 x 10 ⁹ beads/mL
Detector / Sample Diluent	101353	523824	50 nM
Bead Diluent	100458	526507	
SBG Dilution Buffer	100376	515316	
SBG Concentrate	103397	515406	
Bead Conjugation Buffer	100453	524508	
Bead Wash Buffer	101352	521016	
Bead Blocking Buffer	100457	521017	
Biotinylation Reaction Buffer	100460	523109	
EDC, 10mg	100022	RM2025060000020	
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

Dipesh Patel
Name

Sr. QA Specialist
Title

26 SEP 2025
Signature/Date