

Product Number:	102040
Lot Number:	502482
Expiration:	18-Dec-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	015806
Biotinylation Reaction Buffer	101358	015108
Bead Conjugation Buffer	101357	015107
Bead Blocking Buffer	101356	935219
Bead Wash Buffer	101355	015106
Bead Diluent Buffer	100458	007710

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 29 Jul 2020

Name

Title

Signature/Date