

Product Number:	102040
Lot Number:	502470
Expiration:	18-Dec-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	015806
Biotinylation Reaction Buffer	101358	015108
Bead Conjugation Buffer	101357	011706
Bead Blocking Buffer	101356	935219
Bead Wash Buffer	101355	008007
Bead Diluent Buffer	100458	007710

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 20/11/2020
Signature/Date