

Product Number:	102040
Lot Number:	503319
Expiration:	21-Jun-2022

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	134921
Biotinylation Reaction Buffer	101358	117227
Bead Conjugation Buffer	101357	127410
Bead Blocking Buffer	101356	127513
Bead Wash Buffer	101355	127507
Bead Diluent Buffer	100458	122710

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 21Jan2022

Signature/Date