

Product Number:	102040	
Lot Number:	503268	
Expiration:	21-Jun-2022	

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	129126
Biotinylation Reaction Buffer	101358	117227
Bead Conjugation Buffer	101357	127410
Bead Blocking Buffer	101356	127513
Bead Wash Buffer	101355	122510
Bead Diluent Buffer	100458	122710

Review/Approval

*Shivani Goel*

Name

QC

Title

*Shivani Goel 17Dec2021*

Signature/Date