

Product Number:	102040	
Lot Number:	502678	
Expiration:	11-Aug-2021	

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	031906
Biotinylation Reaction Buffer	101358	034007
Bead Conjugation Buffer	101357	034002
Bead Blocking Buffer	101356	024106
Bead Wash Buffer	101355	031113
Bead Diluent Buffer	100458	022412

Review/Approval

Junpil Park

QC Supervisor

 05 Jan 2021

Name

Title

Signature/Date