

Product Number:	102040	
Lot Number:	502690	
Expiration:	28-Aug-2021	

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	031906
Biotinylation Reaction Buffer	101358	025909
Bead Conjugation Buffer	101357	034002
Bead Blocking Buffer	101356	024109
Bead Wash Buffer	101355	031113
Bead Diluent Buffer	100458	035807

Review/Approval

Junpil Park  
Name

QC Supervisor  
Title

 13 Jan 2021  
Signature/Date