

Product Number:	102040
Lot Number:	502248
Expiration:	29-Aug-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	002217
Biotinylation Reaction Buffer	101358	924108
Bead Conjugation Buffer	101357	935108
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	928912
Bead Diluent Buffer	100458	928215

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 20 Feb 2020
Name	Title	Signature/Date