

Product Number:	102040
Lot Number:	502170
Expiration:	19-Jun-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	928214
Biotinylation Reaction Buffer	101358	924104
Bead Conjugation Buffer	101357	920505
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	919213
Bead Diluent Buffer	100458	917022

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	<i>20 Dec 2019</i>
Name	Title	Signature/Date	