

Product Number:	102040
Lot Number:	502223
Expiration:	24-Jul-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	932502
Biotinylation Reaction Buffer	101358	924104
Bead Conjugation Buffer	101357	920505
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	932505
Bead Diluent Buffer	100458	921315

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>p. stadler</i>	03 Feb 2020
<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>	