

Product Number:	102040
Lot Number:	502170
Expiration:	11-Jul-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	929214
Biotinylation Reaction Buffer	101358	924104
Bead Conjugation Buffer	101357	920505
Bead Blocking Buffer	101356	935209
Bead Wash Buffer	101355	919213
Bead Diluent Buffer	100458	921319

Review/Approval

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