

Product Number:	102040
Lot Number:	502049
Expiration:	6-Mar-2020

Component	Part Number	Lot Number
Detector/Sample Diluent	101359	924905
Bead Conjugation Buffer	101357	919109
Bead Diluent Buffer	100458	917020
Biotinylation Reaction Buffer	101358	917620
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	910013

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler 30 Sep 2019</i>
Name	Title	Signature/Date