

<b>Product Number:</b>	102040
<b>Lot Number:</b>	502150
<b>Expiration:</b>	6-Mar-2020

<b>Component</b>	<b>Part Number</b>	<b>Lot Number</b>
Detector/Sample Diluent	101359	926016
Biotinylation Reaction Buffer	101358	924104
Bead Conjugation Buffer	101357	920505
Bead Blocking Buffer	101356	906502
Bead Wash Buffer	101355	919213
Bead Diluent Buffer	100458	917022

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	10 Dec 2019
Name	Title	Signature/Date	