

<b>Product Number:</b>	102040
<b>Lot Number:</b>	503775
<b>Expiration:</b>	5-Dec-2023

<b>Component</b>	<b>Item Number</b>	<b>Lot Number</b>
Detector/Sample Diluent	101359	234016
Biotinylation Reaction Buffer	101358	234206
Bead Conjugation Buffer	101357	233907
Bead Blocking Buffer	101356	234208
Bead Wash Buffer	101355	234302
Bead Diluent Buffer	100458	234304

Review/Approval

Sonia Khillan      QC Manager      Sonia Khillan / 22 DEC 2022  
Name    Title    Signature/Date