

<b>Product Number:</b>	102040
<b>Lot Number:</b>	501680
<b>Expiration:</b>	5-Nov-2019

<b>Component</b>	<b>Item Number</b>	<b>Lot Number</b>
Detector/Sample Diluent	101359	831007
Bead Conjugation Buffer	101357	830909
Bead Diluent Buffer	100458	833308
Biotinylation Reaction Buffer	101358	900922
Bead Blocking Buffer	101356	830905
Bead Wash Buffer	101355	830904

Review/Approval

Patrizia Stadler	QC Assoc Director	<i>P. Stadler</i> 28 Feb 2019
Name	Title	Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.