

Product Number:	101354
Lot Number:	501919
Expiration:	12-Feb-2020

Component	Item Number	Lot Number
Beads	103207	912111
Detector/Sample Diluent	101359	917613
Biotinylation Reaction Buffer	101358	904309
Bead Conjugation Buffer	101357	916503
Bead Blocking Buffer	101356	906514
Bead Wash Buffer	101355	919213
Bead Diluent Buffer	100458	917020

Bead Count (beads/mL)	3.19x10 ⁹
--------------------------	----------------------

Review/Approval

Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i> 18 Jul 2019
_____ Name	_____ Title	_____ Signature/Date

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