

Product Number:	101354
Lot Number:	501606
Expiration:	5-Nov-2019

Component	Item Number	Lot Number
Beads	103207	822906
Detector/Sample Diluent	101359	831007
Biotinylation Reaction Buffer	101358	900922
Bead Conjugation Buffer	101357	830909
Bead Blocking Buffer	101356	830905
Bead Wash Buffer	101355	830904
Bead Diluent Buffer	100458	833308

Bead Count (beads/mL)	3.04x10 ⁹
--------------------------	----------------------

Review/Approval

Patrizia Stadler QC Assoc. Director

p. Stadler 25 Jan 2019

Name

Title

Signature/Date

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