

Product Number:	103217
Lot Number:	820504
Expiration:	24-Jul-2020
Storage:	-80°C

Laboratory Analysis

Concentration (pg/mL)	2565
-----------------------	------

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date

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