

Product Number:	102514
Lot Number:	936010
Expiration:	26-Dec-2021
Storage:	- 20°C

Laboratory Analysis

Concentration (pg/ml)	600
------------------------------	-----

Review/Approval Shwani Goel QC Shwani Goel 22 Oct 2021
Name Title Signature/Date

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