

Product Number:	101735
Lot Number:	501141
Expiration:	19-Dec-2018

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	MF-121917-13
SBG Concentrate	100439	80372

SBG Concentrate Concentration (nM)	119
------------------------------------	-----

Review/Approval

Sivasankari Murugar QC Supervisor
Name Title


Signature/Date 21 FEB 2018

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