

Product Number:	101735
Lot Number:	501218
Expiration:	3-Oct-2018

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	MF-013018-15
SBG Concentrate	100439	72765

SBG Concentrate Concentration (nM)	165
---------------------------------------	-----

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 17 APR 2018
Signature/Date

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