

Product Number:	101735
Lot Number:	501115
Expiration:	28-Nov-2018


Component	Item Number	Lot Number
SBG Dilution Buffer	100376	MF-121917-13
SBG Concentrate	100439	73328

SBG Concentrate Concentration (nM)	128
------------------------------------	-----

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date 07 FEB 2018

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