

Product Number:	102295
Lot Number:	501309
Expiration:	17-Apr-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	81072

SBG Concentrate Concentration (nM)	186
---------------------------------------	-----

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 05/01/2018
Signature/Date

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