

Product Number:	102295
Lot Number:	501750
Expiration:	13-Mar-2020

Component	Item Number	Lot Number
SBG Concentrate	103397	907217

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

01 Apr 2019

Name

Title

Signature/Date

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