

Product Number:	102295
Lot Number:	503337
Expiration:	19-Jul-2022

Component	Item Number	Lot Number
SBG Concentrate	103397	120035

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

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 02 Feb 2022
Signature/Date

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