

Product Number:	103297
Lot Number:	120502
Expiration:	24-Jul-2022
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 26 Jul 2021

Signature/Date