

Product Number:	103297
Lot Number:	116905
Expiration:	18-Jun-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 25 Jun 2021
Signature/Date