

Product Number:	103297
Lot Number:	317909
Control Number:	317909A
Expiration:	28-Jun-2024
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval Sonia Khillan QC Manager Sonia Keller 12JUL2023
Name Title Signature/Date