

Product Number:	103297
Lot Number:	309304
Expiration:	3-Apr-2024
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval Sonia Khilkin QC Manager Sonia Khilkin 06APR2023  
Name Title Signature/Date