

Product Number:	103297
Lot Number:	126103
Expiration:	18-Aug-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 21 Sep 2021
Signature/Date