

<b>Product Number:</b>	103297
<b>Lot Number:</b>	110702
<b>Expiration:</b>	17-Apr-2022
<b>Storage:</b>	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Junpil Park                      QC    Supervisor                      [Signature]                      21 Apr 2021  
Name                                      Title                                      Signature/Date