

Product Number:	103297
Lot Number:	124006
Expiration:	28-Aug-2022
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Junpil Park

QC

Supervisor



01sep2021

Name

Title

Signature/Date