

Product Number:	103297
Lot Number:	523802
Expiration:	26-Feb-2027
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-X	PASS

Review/Approval

Hardi Caswell
Name

Director, QC
Title

Hardi Caswell 08-Sep-2025
Signature/Date