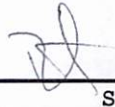


Product Number:	103297
Lot Number:	507608
Expiration:	17-Sep-2026
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval	Robert Somogyi	QC Manager			2024.09.05
	Name	Title	Signature/Date		