

Product Number:	103297
Lot Number:	422201
Expiration:	9-Feb-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

 16 AUG 2024

Name

Title

Signature/Date