


Product Number:	103297
Lot Number:	408502
Expiration:	25-Sep-2025
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval ROBERT SOMOGYI QC MANAGER  01 APR 2024

Name	Title	Signature/Date
------	-------	----------------