

Product Number:	103397
Lot Number:	401702
Expiration:	17-Jul-2025
Storage:	2-8°C

Laboratory Analysis		
Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD(X/1) Analyzer	PASS

RESULT	
SBG Concentrate Concentration (nM)	50

Review/Approval

ROBERT SOMOGYI

QC MANAGER

 31 JAN 2024

Name

Title

Signature/Date