

Product Number:	103397
Lot Number:	319211
Control Lot Number:	319211A
Expiration:	8-Sep-2025
Storage:	2-8°C

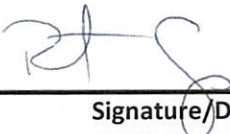
Laboratory Analysis		
Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD(X/1) Analyzer	PASS

RESULT	
SBG Concentrate Concentration (nM)	50

Review/Approval

ROBERT SOMOGYI

QC MANAGER

12 MAR 2024

Name

Title

Signature/Date