

Product Number:	103397
Lot Number:	319213
Control Lot Number:	319213A
Expiration:	4-Mar-2026
Storage:	2-8°C

Laboratory Analysis

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

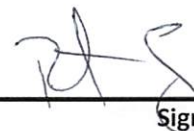
RESULT

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval

Robert Somogyi

QC Manager



09 SEP 2024

Name

Title

Signature/Date