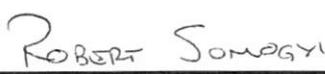
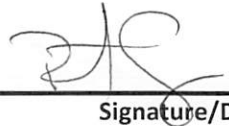


Product Number:	104433
Lot Number:	319214
Control Lot Number:	319214A
Expiration:	11-Jul-2026
Storage:	-20°C

Laboratory Analysis		
Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD-X Analyzer	PASS

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

Review/Approval

Name: ROBERT SOMOGYI Title: QC MANAGER Signature/Date: [Signature] 11 SEP 2024