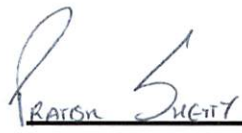


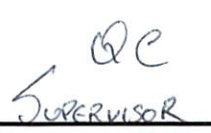
Product Number:	102260
Lot Number:	211107
Expiration:	21-Apr-2024
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	(ug/mL)
HD-X Analyzer	Concentration	437

Review/Approval


RATON SUGIT
Name


QC SUPERVISOR
Title


20 APR 2023
Signature/Date