

Product Number:	102260
Lot Number:	312601
Control Number:	312601A
Expiration:	6-May-2025
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-X Analyzer	Concentration	290

Review/Approval





Name: Pratik Smetty Title: Q. Supervisor Signature/Date: Pratik Smetty 22 MAY 2023