

Product Number:	103710
Lot Number:	415013
Expiration:	29-May-2026
Storage:	-80°C
Total Antigen Concentration:	1.00 ug/mL
Phosphorylated Antigen Concentration:	0.70 ug/mL

Laboratory Analysis

Means of Analysis	Characteristic	(ug/mL)
HD-X Analyzer	Concentration	0.70

Review/Approval

Robert Somogyi
Name

QC Manager
Title

 18 FEB 2025
Signature/Date