

Product Number:	18910024
Lot Number:	513654000002540
Expiration Date:	19Apr2029
Storage:	-20°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Peptide Molecular Weight	Mass Spectral Analysis	PASS

Review/Approval

Dipesh Patel

Sr. Doc Control - QA

at 18 Jun 2025

Name

Title

Signature/Date