

Means of Analysis	PASS / FAIL
HD-X	PASS



DP 28 Mar 2025
Signature/Date