



## CERTIFICATE OF ANALYSIS

**Product Description:** Human TSLP Calibrator  
**Catalog Number:** C01D3-2  
**Lot Number:** A00S0120  
**Expiration Date:** 31 OCT 2029  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
TSLP	580000	fg/mL

**Reconstitution Volume:** 1000 µL

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		13 NOV 2024