



## CERTIFICATE OF ANALYSIS

**Product Description:** TH17 Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0085-2  
**Lot Number:** A0080366  
**Expiration Date:** 31 MAY 2028  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A	3230	pg/mL
IL-21	1020	pg/mL
IL-31	1010	pg/mL
IL-27	15900	pg/mL
IL-23	5270	pg/mL
IL-22	521	pg/mL
MIP-3 $\alpha$	492	pg/mL

**Reconstitution Volume:** 1000  $\mu$ L  $\mu$ L

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		11 JUL 2023