



CERTIFICATE OF ANALYSIS

Product Description: Cytokine Panel 2 (human) Calibrator Blend
Catalog Number: C0084-2
Lot Number: A0080444
Expiration Date: 30 SEP 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A/F	6150	pg/mL
IL-17B	1230	pg/mL
IL-17C	2770	pg/mL
IL-17D	9010	pg/mL
IL-1RA	980	pg/mL
IL-3	3510	pg/mL
IL-9	1550	pg/mL
TSLP	521	pg/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
Review/Approval	Karen Hamilla	Quality		05 NOV 2025