

**CERTIFICATE OF ANALYSIS**

Product Description: Cytokine Panel 2 (human) Calibrator Blend
Catalog Number: C0084-2
Lot Number: A0080346
Expiration Date: 31 DEC 2027
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A/F	6000	pg/mL
IL-17B	1390	pg/mL
IL-17C	2510	pg/mL
IL-17D	8210	pg/mL
IL-1RA	1040	pg/mL
IL-3	3010	pg/mL
IL-9	1560	pg/mL
TSLP	475	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	Nithya Santhanam	Quality		10 FEB 2023