



CERTIFICATE OF ANALYSIS

Product Description: TH17 Panel 1 (human) Calibrator Blend
Catalog Number: C0085-2
Lot Number: A0080354
Expiration Date: 31 DEC 2027
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A	2850	pg/mL
IL-21	1030	pg/mL
IL-31	1070	pg/mL
IL-27	17800	pg/mL
IL-23	5160	pg/mL
IL-22	494	pg/mL
MIP-3 α	495	pg/mL

Reconstitution Volume: 1000 μ 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
Review/Approval	Karen Hamilla	Quality		03 MAR 2023