



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: TH17 Panel 1 (human) Calibrator Blend

Catalog Number: C0085-2

Lot Number: A0080263

Expiration Date: 30 SEP 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A	3110	pg/mL
IL-21	932	pg/mL
IL-31	970	pg/mL
IL-27	19900	pg/mL
IL-23	5110	pg/mL
IL-22	521	pg/mL
MIP-3 α	506	pg/mL

Reconstitution Volume: 1000 μ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 FEB 2021