



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** TH17 Panel 1 (human) Calibrator Blend

**Catalog Number:** C0085-2

**Lot Number:** A0080241

**Expiration Date:** 28 FEB 2025

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A	3140	pg/mL
IL-21	1180	pg/mL
IL-31	964	pg/mL
IL-27	18200	pg/mL
IL-23	5540	pg/mL
IL-22	542	pg/mL
MIP-3 $\alpha$	508	pg/mL

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

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		10 MAR 2020