



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0085-2

**Lot Number:** A0080232

**Expiration Date:** 31 OCT 2024

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A	2570	pg/mL
IL-31	1050	pg/mL
IL-27	21000	pg/mL
IL-23	5730	pg/mL
IL-22	527	pg/mL
MIP-3 $\alpha$	565	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
Review/Approval	Karen Hamilla	Quality		15 NOV 2019