



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0246-2

**Lot Number:** A0080192

**Expiration Date:** 31 MAY 2024

**Storage Temperature:** 2-8°C

### Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
MIP-3 $\alpha$	1 150	pg/mL
IL-22	544	pg/mL
IL-23	8 330	pg/mL
IL-17C	1 720	pg/mL
IL-31	33 400	pg/mL
IL-17F	88 500	pg/mL
IL-16	2 000	pg/mL
IL-17A	566	pg/mL
IL-17E/IL-25	4 600	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		27 SEP 2019