



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: TH17 Panel 1 (human) Calibrator Blend

Catalog Number: C0085-2

Lot Number: A0080219

Expiration Date: 30 APR 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A	2 640	pg/mL
IL-21	1 200	pg/mL
IL-31	1 060	pg/mL
IL-27	19 300	pg/mL
IL-23	5 220	pg/mL
IL-22	545	pg/mL
MIP-3a	393	pg/mL

Reconstitution Volume: 1000 µL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 JUN 2019