



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: TH17 Panel 1 (human) Calibrator Blend

Catalog Number: C0085-2

Lot Number: A0080209

Expiration Date: 31 MAY 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A	2720	pg/mL
IL-21	1370	pg/mL
IL-31	1040	pg/mL
IL-27	12100	pg/mL
IL-23	4830	pg/mL
IL-22	533	pg/mL
MIP-3 α	587	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	Dawn Stollar	Quality		29 MAY 2019