



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: TH17 Panel 1 (mouse) Calibrator Blend

Catalog Number: C0246-2

Lot Number: A0080220

Expiration Date: 30 JUN 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
MIP-3 α	988	pg/mL
IL-22	488	pg/mL
IL-23	8420	pg/mL
IL-17C	1620	pg/mL
IL-31	30000	pg/mL
IL-21	10400	pg/mL
IL-17F	84100	pg/mL
IL-16	2000	pg/mL
IL-17A	559	pg/mL
IL-17E/IL-25	4840	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	Karen Hamilla	Quality		24 JUL 2019