



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: TH17 Panel 1 (mouse) Calibrator Blend

Catalog Number: C0246-2

Lot Number: A0080146

Expiration Date: 31 AUG 2022

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	ms MIP-3 α	ms IL-22	ms IL-23	ms IL-17C	ms IL-31	ms IL-21	ms IL-17F	ms IL-16	ms IL-17A	ms IL-17E/IL-25
Reconstituted Calibrator (pg/mL)	1130	510	8380	1770	28 500	12 000	85 000	2420	550	4540

Reconstitution Volume: 1000 μ L

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 SEP 2017