



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0246-2

**Lot Number:** A0080154

**Expiration Date:** 31 OCT 2022

**Storage Temperature:** 2-8°C

**Concentration of Materials Tested:**

	Assay									
	ms MIP-3α	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
<b>Reconstituted Calibrator (pg/mL)</b>	1060	555	8690	1500	30 700	16 100	87 800	2030	585	5780

**Reconstitution Volume:** 1000 µL

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	27 OCT 2017