



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: TH17 Panel 1 (human) Calibrator Blend

Catalog Number: C0085-2

Lot Number: A0080206

Expiration Date: 31 JAN 2022

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay						
	hu IL-17A	hu IL-21	hu IL-31	hu IL-27	hu IL-23	hu IL-22	hu MIP-3α
Reconstituted Calibrator (pg/mL)	2900	842	962	16 100	4600	551	566

Reconstitution Volume: 1000 µL

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		28 MAR 2019