



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07515

Product Description: V-PLEX TH17 Panel 1 (human) Calibrator Blend
Catalog Number: C0085-2
Lot Number: A0080101
Expiration Date: 30 SEP 2021
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)	2730	1040	881	17100	4900	560	449

Reconstitution Volume: 1000 µL
Reconstituted Concentration: 1X

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. 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>	31 JAN 2017