



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07643

Product Description: V-PLEX TH17 Panel 1 (human) Calibrator Blend
Catalog Number: C0085-2
Lot Number: A0080111
Expiration Date: 30 NOV 2021
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay						
	hu IL-17A/F	hu IL-21	hu IL-31	hu IL-27	hu IL-23	hu IL-22	hu MIP-3α
Reconstituted Calibrator (pg/mL)	2600	913	895	15100	4350	523	443

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>	16 DEC 2016