



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06306

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080069

Expiration Date: 31 OCT 2020

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	hu GM-CSF	hu IL-1 α	hu IL-5	hu IL-7	hu IL-12/23 p40	hu IL-15	hu IL-16	hu IL-17A	hu TNF- β	hu VEGF
Reconstituted Calibrator (pg/mL)	996	405	912	838	3220	779	2530	6010	632	1020

Reconstitution Volume: 1.0 mL

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>	14DEC2015