



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05351

Product Description: Cytokine Panel 1 (human) Calibrator Blend
Catalog Number: C0050-2
Lot Number: A0080051
Expiration Date: 29 FEB 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)	963	350	892	650	2840	654	2320	5500	580	941

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	Dana Short	Quality Control		02 APR 2015