



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06120

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080068

Expiration Date: 30 SEP 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)	972	388	891	836	3110	762	2580	5890	625	1070

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	Monika Jain	Quality Assurance	<i>Monika Jain</i>	28 OCT 2015