



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05075

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080047

Expiration Date: 31 DEC 2019

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)	978	368	883	640	2840	673	2450	5940	591	961

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	Caitlin Hansen	Quality Control		13 JAN 2015