



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04572

**Product Description:** Cytokine Panel 1 (human) Calibrator Blend

**Catalog Number:** C0050-2

**Lot Number:** A0080041

**Expiration Date:** 30 JUN 2019

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay									
	hu GM-CSF	hu IL-1 $\alpha$	hu IL-5	hu IL-7	hu IL-12/23 p40	hu IL-15	hu IL-16	hu IL-17A	hu TNF- $\beta$	hu VEGF
Reconstituted Calibrator (pg/mL)	950	358	817	628	3070	683	2810	5670	581	996

**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
<b>Review/Approval</b>	Caitlin Hansen	Quality Control		23 JUL 2014