



### CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05185

**Product Description:** Cytokine Panel 1 (human) Control Pack  
**Catalog Number:** C4050-1  
**Lot Number:** CP000026  
**Expiration Date:** 31 OCT 2019

**Concentration of Materials Tested:**

Units = pg/mL	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 Control 1 (A00C0280)	346	252	350	315	1310	191	1340	2020	221	254
Reconstituted Control 2 (A00C0281)	53.9	55.8	113	46.7	202	37	366	281	34.4	92.4
Reconstituted Control 3 (A00C0300)	7.62	17.3	36.3	6.81	30.1	6.63	119	50.1	4.94	32.6

**Reconstitution Volume:** 250  $\mu$ L

Control concentrations were assigned across three kit lots.

Cytokine Panel 1 (human) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0280	2-8 $^{\circ}$ C	31 OCT 2019
Cytokine Panel 1 (human) Control 2	A00C0281	2-8 $^{\circ}$ C	31 OCT 2019
Cytokine Panel 1 (human) Control 3	A00C0300	2-8 $^{\circ}$ C	31 JAN 2020

**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>	Dana Short	Quality Control		05 FEB 2015