



# CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05707

**Product Description:** Cytokine Panel 1 (human) Control Pack

**Catalog Number:** C4050-1

**Lot Number:** CP000030

**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 (A00C0325)	362	167	371	241	1290	241	1160	2590	196	373
Reconstituted Control 2 (A00C0326)	55.8	57.8	123	42.0	208	36.5	328	327	31.7	127
Reconstituted Control 3 (A00C0282)	8.84	18.2	38.3	7.09	31.1	6.97	113	37.7	4.79	21.3

**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	A00C0325	2-8°C	30 APR 2020
Cytokine Panel 1 (human) Control 2	A00C0326	2-8°C	31 MAY 2020
Cytokine Panel 1 (human) Control 3	A00C0282	2-8°C	31 OCT 2019

**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>	Anastasia Belilovets	Quality Assurance		10 Jul 2015