



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4050-1

**Lot Number:** CP000065

**Expiration Date:** 31 OCT 2022

### 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
Control 1 (A00C0528)	374	178	322	348	1380	288	1210	2350	229	428
Control 2 (A00C0529)	64.7	62.3	105	56.4	230	46.3	329	318	41.0	135
Control 3 (A00C0530)	9.35	18.3	33.7	8.49	34.7	8.87	104	47.7	5.97	45.1

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

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0528	2-8°C	31 OCT 2022
Cytokine Panel 1 (human) Control 2	A00C0529	2-8°C	31 OCT 2022
Cytokine Panel 1 (human) Control 3	A00C0530	2-8°C	31 OCT 2022

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	16 AUG 2018