



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (human) Control Pack

Catalog Number: C4050-1

Lot Number: CP000080

Expiration Date: 31 AUG 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

Units = pg/mL	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
Control 1 (A00C0610)	408	183	343	346	1430	286	1200	2020	247	384
Control 2 (A00C0617)	64.1	58.1	109	58.0	204	45.0	334	277	37.9	114
Control 3 (A00C0612)	9.70	16.3	32.7	7.79	32.6	8.00	100	42.9	5.75	37.5

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0610	2-8°C	31 AUG 2023
Cytokine Panel 1 (human) Control 2	A00C0617	2-8°C	30 SEP 2023
Cytokine Panel 1 (human) Control 3	A00C0612	2-8°C	31 AUG 2023

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	08 NOV 2018