



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4050-1

**Lot Number:** CP000050

**Expiration Date:** 31 DEC 2021

### 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 (A00C0454)	388	203	343	386	1580	302	1420	2640	261	429
Control 2 (A00C0455)	60.6	61.9	108	59.2	236	53.0	363	387	43.8	136
Control 3 (A00C0456)	9.33	20.3	32.8	9.41	39.4	9.62	106	54.4	5.54	49.0

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

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 1 (human) Control 1	A00C0454	2-8°C	31 DEC 2021
Cytokine Panel 1 (human) Control 2	A00C0455	2-8°C	31 DEC 2021
Cytokine Panel 1 (human) Control 3	A00C0456	2-8°C	31 DEC 2021

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>	25 APR 2017