



### CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4049-1

**Lot Number:** CP000069

**Expiration Date:** 28 FEB 2023

#### Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu IFN- $\gamma$	hu IL-1 $\beta$	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70	hu IL-13	hu TNF- $\alpha$
Reconstituted Control 1 (A00C0531)	582	208	407	101	271	233	125	194	168	89.5
Reconstituted Control 2 (A00C0532)	164	52.7	42.1	20.6	43.5	37.4	24.2	36.4	48.8	17.4
Reconstituted Control 3 (A00C0533)	40.9	10.6	3.54	2.95	8.03	6.13	3.49	7.29	13.4	2.21

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

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0531	2-8°C	31 MAR 2023
Proinflammatory Panel 1 (human) Control 2	A00C0532	2-8°C	28 FEB 2023
Proinflammatory Panel 1 (human) Control 3	A00C0533	2-8°C	31 MAR 2023

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>	10 SEP 2018