



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Control Pack

Catalog Number: C4049-1

Lot Number: CP000079

Expiration Date: 31 AUG 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

Units = pg/mL	Assay									
	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70	hu IL-13	hu TNF- α
Reconstituted Control 1 (A00C0614)	512	203	390	102	264	235	133	186	170	92.0
Reconstituted Control 2 (A00C0615)	132	47.1	38.3	20.0	46.8	36.2	25.3	35.6	43.8	17.2
Reconstituted Control 3 (A00C0616)	40.9	10.9	3.34	3.05	8.18	5.93	3.32	6.99	12.7	2.53

Reconstitution Volume: 250 μ L

Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0614	2-8°C	31 AUG 2023
Proinflammatory Panel 1 (human) Control 2	A00C0615	2-8°C	31 AUG 2023
Proinflammatory Panel 1 (human) Control 3	A00C0616	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>	21 SEP 2018