



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06177

Product Description: Proinflammatory Panel 1 (human) Control Pack

Catalog Number: C4049-1

Lot Number: CP000034

Expiration Date: 31 AUG 2019

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 (A00C0347)	535	199	390	70.3	276	180	174	143	209	54.0
Reconstituted Control 2 (A00C0343)	140	47.1	37.0	15.0	53.6	28.3	26.1	28.0	54.6	12.4
Reconstituted Control 3 (A00C0275)	91.5	15.1	18.4	3.41	22.9	10.3	7.67	7.32	44.4	3.50

Reconstitution Volume: 250 μ L

Control concentrations were assigned across three kit lots.

Proinflammatory Panel 1 (human) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0347	2-8°C	31 MAY 2020
Proinflammatory Panel 1 (human) Control 2	A00C0343	2-8°C	31 MAY 2020
Proinflammatory Panel 1 (human) Control 3	A00C0275	2-8°C	31 AUG 2019

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Tanisha Lucas	Quality Assurance	<i>Tanisha Lucas</i>	05 NOV 2015