



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06159

Product Description: Proinflammatory Panel 1 (rat) Control Pack
Catalog Number: C4044-1
Lot Number: CP000035
Expiration Date: 31 OCT 2020

Concentration of Materials Tested:

Units = pg/mL	Assay									
	rt IFN- γ	rt IL-1 β	rt IL-2	rt IL-4	rt IL-5	rt IL-6	rt KC/GRO	rt IL-10	rt IL-13	rt TNF- α
Reconstituted Control 1 (A00C0348)	5930	12400	N/A	951	3380	8090	554	19800	1580	803
Reconstituted Control 2 (A00C0349)	1390	7940	N/A	312	1730	2880	327	4360	495	282
Reconstituted Control 3 (A00C0350)	736	1820	N/A	81.5	770	726	158	1640	162	86.9

Reconstitution Volume: 250 μ L

Control concentrations were assigned across three kit lots.

Proinflammatory Panel 1 (rat) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0348	2-8°C	31 OCT 2020
Proinflammatory Panel 1 (rat) Control 2	A00C0349	2-8°C	31 OCT 2020
Proinflammatory Panel 1 (rat) Control 3	A00C0350	2-8°C	31 OCT 2020

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	Monika Jain	Quality Assurance	<i>Monika Jain</i>	02 NOV 2015