



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04627

**Product Description:** Proinflammatory Panel 1 (rat) Control Pack**Catalog Number:** C4044-1**Lot Number:** CP000019**Expiration Date:** 31 JUL 2019**Concentration of Materials Tested:**

Units = pg/mL	Assay									
	rt IFN- $\gamma$	rt IL-1 $\beta$	rt IL-2	rt IL-4	rt IL-5	rt IL-6	rt KC/GRO	rt IL-10	rt IL-13	rt TNF- $\alpha$
Reconstituted Control 1 (A00C0257)	8510	34400	N/A	1950	7660	16500	719	41500	2400	1440
Reconstituted Control 2 (A00C0258)	1550	11800	N/A	460	2560	3290	535	6240	691	457
Reconstituted Control 3 (A00C0259)	628	2340	N/A	112	831	842	193	2080	177	164

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

Control concentrations were assigned across 3 kit lots.

Proinflammatory Panel 1 (rat) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (rat) Control 1	A00C0257	2-8°C	31 JUL 2019
Proinflammatory Panel 1 (rat) Control 2	A00C0258	2-8°C	31 JUL 2019
Proinflammatory Panel 1 (rat) Control 3	A00C0259	2-8°C	31 JUL 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
<b>Review/Approval</b>	Monika Jain	Quality Assurance	<i>Monika Jain</i>	21 AUG 2015