



# CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06866

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

**Catalog Number:** C4049-1

**Lot Number:** CP000038

**Expiration Date:** 28 FEB 2021

## 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 (A00C0384)	647	207	418	99.6	265	242	131	161	149	95.5
Reconstituted Control 2 (A00C0385)	183	49.5	33.6	18.1	54.5	39.3	26.7	31.1	43.3	16.4
Reconstituted Control 3 (A00C0386)	47.8	9.90	3.40	5.02	7.72	6.78	3.45	5.56	12.7	2.17

**Reconstitution Volume:** 250  $\mu$ 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	A00C0384	2-8°C	28 FEB 2021
Proinflammatory Panel 1 (human) Control 2	A00C0385	2-8°C	31 MAR 2021
Proinflammatory Panel 1 (human) Control 3	A00C0386	2-8°C	31 MAR 2021

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>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	06 MAY 2016