



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C4049-1

**Lot Number:** CP000068

**Expiration Date:** 31 DEC 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 (A00C0448)	682	215	431	101	275	223	133	168	150	89.6
Reconstituted Control 2 (A00C0532)	164	52.7	42.1	20.6	43.5	37.4	24.2	36.4	48.8	17.4
Reconstituted Control 3 (A00C0450)	52.3	11	3.83	2.88	9.03	5.91	3.61	6.12	12.3	2.72

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

Control concentrations were assigned across three kit lots.

Proinflammatory Panel 1 (mouse) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Control 1	A00C0448	2-8° C	31 DEC 2021
Proinflammatory Panel 1 (human) Control 2	A00C0532	2-8° C	28 FEB 2023
Proinflammatory Panel 1 (human) Control 3	A00C0450	2-8° C	31 DEC 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>	10 SEP 2018