



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07933

Product Description: Proinflammatory Panel 1 (human) Control Pack
Catalog Number: C4049-1
Lot Number: CP000045
Expiration Date: 31 DEC 2021

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 (A00C0448)	682	215	431	101	275	223	133	168	150	89.6
Reconstituted Control 2 (A00C0449)	183	52.0	42.3	20.3	50.7	40.4	27.8	33.2	46.7	18.0
Reconstituted Control 3 (A00C0450)	52.3	11.0	3.83	2.88	9.03	5.91	3.61	6.12	12.3	2.72

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	A00C0448	2-8°C	31 DEC 2021
Proinflammatory Panel 1 (human) Control 2	A00C0449	2-8°C	31 JAN 2022
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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 FEB 2017