



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05050

Product Description: Proinflammatory Panel 1 (mouse) Control Pack

Catalog Number: C4048-1

Lot Number: CP000025

Expiration Date: 31 OCT 2019

Concentration of Materials Tested:

Units = pg/mL	Assay									
	ms IFN- γ	ms IL-1 β	ms IL-2	ms IL-4	ms IL-5	ms IL-6	ms KC/GRO	ms IL-10	ms IL-12p70	ms TNF- α
Reconstituted Control 1 (A00C0289)	479	623	933	505	381	1740	902	1140	13100	172
Reconstituted Control 2 (A00C0290)	26.8	48.7	67.8	89.3	46.0	271	115	283	3060	32.5
Reconstituted Control 3 (A00C0291)	3.05	1.57	7.43	24.6	7.83	66.0	15.3	115	1010	5.19

Reconstitution Volume: 250 μ 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 (mouse) Control 1	A00C0289	2-8° C	31 OCT 2019
Proinflammatory Panel 1 (mouse) Control 2	A00C0290	2-8° C	31 DEC 2019
Proinflammatory Panel 1 (mouse) Control 3	A00C0291	2-8° C	31 OCT 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
Review/Approval	Caitlin Hansen	Quality Control		19 DEC 2014