



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05836

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend

Catalog Number: C0048-2

Lot Number: A0080060

Expiration Date: 31 JUL 2020

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	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 Calibrator (pg/mL)	878	1460	2310	1540	850	5200	1880	2630	31 600	536

Reconstitution Volume: 1.0 mL

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	Monika Jain	Quality Assurance	<i>Monika Jain</i>	03Aug2015