



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06000

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080061

Expiration Date: 30 JUN 2020

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	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 Calibrator (pg/mL)	1450	478	1350	N/A	615	502	327	431	489	300

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>	22 SEP 2015