



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04622

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080040

Expiration Date: 31 MAY 2019

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)	1460	495	1460	198	767	495	334	421	466	312

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	Caitlin Hansen	Quality Control		11 AUG 2014