



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06131

**Product Description:** Proinflammatory Panel 1 (human) Calibrator Blend

**Catalog Number:** C0049-2

**Lot Number:** A0080070

**Expiration Date:** 30 SEP 2020

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay									
	hu IFN- $\gamma$	hu IL-1 $\beta$	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10	hu IL-12p70	hu IL-13	hu TNF- $\alpha$
Reconstituted Calibrator (pg/mL)	1500	496	1380	262	653	575	312	425	539	327

**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
<b>Review/Approval</b>	Monika Jain	Quality Assurance	<i>Monika Jain</i>	28 OCT 2015