



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05231

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

**Catalog Number:** C0049-2

**Lot Number:** A0080049

**Expiration Date:** 31 JAN 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)	1340	515	1290	183	650	505	301	397	490	295

**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>	Dana Short	Quality Control		20 FEB 2015