



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05051

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

**Catalog Number:** C0049-2

**Lot Number:** A0080045

**Expiration Date:** 30 NOV 2019

**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)	1510	520	1460	207	749	546	313	404	489	311

**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>	Caitlin Hansen	Quality Control		19 DEC 2014