



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06778

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

**Catalog Number:** C0049-2

**Lot Number:** A0080076

**Expiration Date:** 30 APR 2021

**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)	1650	563	1550	229	751	626	355	480	502	352

**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>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	12 APR 2016