



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05562

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

**Catalog Number:** C0044-2

**Lot Number:** A0080059

**Expiration Date:** 31 MAY 2020


**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay									
	rt IFN- $\gamma$	rt IL-1 $\beta$	rt IL-2	rt IL-4	rt IL-5	rt IL-6	rt KC/GRO	rt IL-10	rt IL-13	rt TNF- $\alpha$
Reconstituted Calibrator (pg/mL)	4820	10 700	N/A	1070	12 100	15 300	2960	26 800	1800	1390

**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>	Anastasia Belilovets	Quality Assurance		16 Jun 2015