



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07207

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

**Catalog Number:** C0044-2

**Lot Number:** A0080083

**Expiration Date:** 31 MAY 2021

**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)	6300	13 700	N/A	1310	9280	12 200	4060	26 100	1670	1340

**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>	11 AUG 2016