



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06475

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

**Catalog Number:** C0044-2

**Lot Number:** A0080066

**Expiration Date:** 31 JAN 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)	5440	13 000	N/A	1280	10 700	13 500	2920	29 000	1810	1440

**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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	08FEB2016