



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06685

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

**Catalog Number:** C0048-2

**Lot Number:** A0080075

**Expiration Date:** 28 FEB 2021

**Storage Temperature:** 2-8°C

### Concentration of Materials Tested:

	Assay									
	ms IFN- $\gamma$	ms IL-1 $\beta$	ms IL-2	ms IL-4	ms IL-5	ms IL-6	ms KC/GRO	ms IL-10	ms IL-12p70	ms TNF- $\alpha$
Reconstituted Calibrator (pg/mL)	938	1670	2630	1660	967	5720	1980	3410	32 200	627

**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>	24 MAR 2016