



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06684

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend
Catalog Number: C0044-2
Lot Number: A0080072
Expiration Date: 28 FEB 2021
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	rt IFN- γ	rt IL-1 β	rt IL-2	rt IL-4	rt IL-5	rt IL-6	rt KC/GRO	rt IL-10	rt IL-13	rt TNF- α
Reconstituted Calibrator (pg/mL)	6550	14 100	N/A	1350	9270	12 300	3650	29 300	1760	1300

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>	24 MAR 2016