



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05511

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend

Catalog Number: C0044-2

Lot Number: A0080056

Expiration Date: 30 APR 2020

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)	6650	N/A	N/A	1070	11 200	13 400	3030	N/A	1850	1380

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	Anastasia Belilovets	Quality Assurance		26 May 2015