



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04160

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend

Catalog Number: C0044-2

Lot Number: A0080035

Expiration Date: 31Oct2018

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)	4890	12 100	N/A	932	10 800	11 800	2800	18 600	1470	1060

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	Caitlin Hansen	Quality Control		19Feb2014