



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05355

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080053
Expiration Date: 31 JUL 2019
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	ms IFN- γ	ms IL-1 β	ms IL-2	ms IL-4	ms IL-5	ms IL-6	ms KC/GRO	ms IL-10	ms IL-12p70	ms TNF- α
Reconstituted Calibrator (pg/mL)	890	1500	2180	1500	869	4200	1840	2800	29 600	532

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	Dana Short	Quality Control		03 APR 2015