



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07646

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080110
Expiration Date: 31 DEC 2021
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)	933	1720	2720	1700	961	5500	1920	3070	31 500	653

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>	16 DEC 2016