



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend

Catalog Number: C0048-2

Lot Number: A0080168

Expiration Date: 31 MAY 2023

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
Reconstitution Calibrator (pg/mL)	933	1430	2420	1610	903	4730	1830	3480	32 600	629

Reconstitution Volume: 1000 μ L

Statement: 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 MAY 2018