



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend

Catalog Number: C0048-2

Lot Number: A0080152

Expiration Date: 30 NOV 2022

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)	644	1510	2980	1670	947	4990	1870	3530	30 400	662

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>	26 JAN 2018