



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080178

Expiration Date: 31 AUG 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	hu IFN- γ	hu IL-1 β	hu IL-2	hu IL-4	hu IL-6	hu IL-8	hu IL-10	hu IL- 12p70	hu IL-13	hu TNF- α
Reconstituted Calibrator (pg/mL)	1430	576	1470	243	764	621	340	542	525	369

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>	11 SEP 2018