



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080144

Expiration Date: 31 AUG 2022

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	hu GM-CSF	hu IL-1 α	hu IL-5	hu IL-7	hu IL-12/23p40	hu IL-15	hu IL-16	hu IL-17A	hu IL-TNF- β	hu VEGF
Reconstituted Calibrator (pg/mL)	1100	459	877	881	3690	842	2800	6480	733	1490

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>	27 SEP 2017