



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080227

Expiration Date: 31 OCT 2024

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
GM-CSF	808	pg/mL
IL-1 α	457	pg/mL
IL-5	895	pg/mL
IL-7	819	pg/mL
IL-12/IL-23p40	3900	pg/mL
IL-15	976	pg/mL
IL-16	3040	pg/mL
IL-17A	5770	pg/mL
TNF- β	654	pg/mL
VEGF	1770	pg/mL

Reconstitution Volume: 1000 μ L

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Carissa Hubbard	Quality		26 DEC 2019