



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 2 (human) Calibrator Blend

Catalog Number: C0084-2

Lot Number: A0080251

Expiration Date: 30 APR 2025

Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-17A/F	6590	pg/mL
IL-17B	1790	pg/mL
IL-17C	2800	pg/mL
IL-17D	9050	pg/mL
IL-1RA	1240	pg/mL
IL-3	3710	pg/mL
IL-9	1620	pg/mL
TSLP	499	pg/mL

Reconstitution Volume: 1000 µL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 MAY 2020