



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 2 (human) Calibrator Blend

Catalog Number: C0084-2

Lot Number: A0080114

Expiration Date: 31 DEC 2021

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay							
	hu IL-17A/F	hu IL-17B	hu IL-17C	hu IL-17D	hu IL-1RA	hu IL-3	hu IL-9	hu TSLP
Reconstituted Calibrator (pg/mL)	5430	1240	2510	7890	1150	3360	1490	555

Reconstitution Volume: 1.0 mL

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>	28 JUN 2017