



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07824

Product Description: V-PLEX Chemokine Panel 1 (human) Calibrator Blend

Catalog Number: C0047-2

Lot Number: A0080080

Expiration Date: 31 MAY 2021

Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	hu Eotaxin	hu MIP-1 β	hu Eotaxin-3	hu TARC	hu IP-10	hu MIP-1 α	hu IL-8	hu MCP-1	hu MDC	hu MCP-4
Reconstituted Calibrator (pg/mL)	1770	1150	6970	1790	2580	1050	83000	497	11600	730

Reconstitution Volume: 1000 μ L

Reconstituted Concentration: 1X

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. 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 JAN 2017