



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 20

Catalog Number: C0328-2

Lot Number: A00U0160

Expiration Date: 30 JUN 2025

Storage Temperature: 2-8°C

Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
CD276/B7-H3	99250	pg	397000	pg/mL
CD20	182000	pg	728000	pg/mL
CD27	8450	pg	33800	pg/mL
CD28	320000	pg	1280000	pg/mL
CTLA-4	3375	pg	13500	pg/mL
GITRL/TNFSF18	2600	pg	10400	pg/mL
OX40/TNFRSF4	4150	pg	16600	pg/mL
PD1 (epitope 1)	2500	pg	10000	pg/mL
PD1 (epitope 2)	2500	pg	10000	pg/mL
TIGIT	7925	pg	31700	pg/mL
TLR1	90000	pg	360000	pg/mL

Reconstitution Volume: 250 µL

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 JUN 2020