



CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 32  
**Catalog Number:** C0719-2  
**Lot Number:** A00U0348  
**Expiration Date:** 31 OCT 2028  
**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
BTLA	230000	pg	920000	pg/mL
CD80/B7-1	16725	pg	66900	pg/mL
IDO-1	42000	pg	168000	pg/mL
proMMP-10	5925	pg	23700	pg/mL
MMP-10 (total)	5925	pg	23700	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
<b>Review/Approval</b>	Karen Hamilla	Quality		06 Feb 2024