



ImmunoSmol

Arginine ELISA Kit

IS-I-0400

Our Arginine (ARG) ELISA kit allows for the determination of ARG in serum and plasma samples in a minimal sample volume of 20µL. The kit is easy to implement and well suited for both preclinical and clinical studies.

SCIENTIFIC BACKGROUND

Arginine (ARG) is a basic non-essential amino acid well known to participate in physiological functions such as **i)** vascular tone homeostasis as contributing to the production of nitric oxide (NO) and **ii)** immune response control, through **NO synthase** (NOS) and **Arginase** enzymes, respectively.

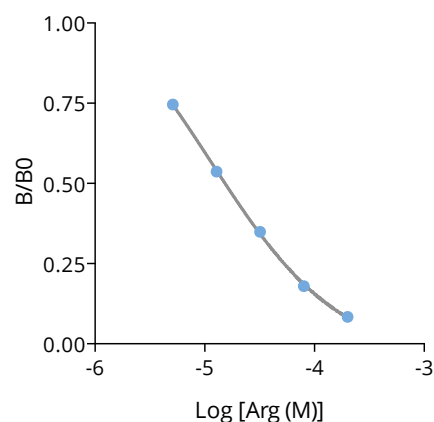
A dysregulation in ARG catabolism has been already linked to tumor progression and immune escape. Indeed, overexpression of Arginase in tumor and immunosuppressive cells (eg. **Myeloid Derived Suppressor Cells**) contributes to an excessive use of ARG thereby limiting its availability for effector immune cells. Arginase represents therefore an attractive therapeutic target to restore **anti-tumor immunity** and inhibitors (eg. INCB001158) are currently under clinical development in several cancer indications. ARG measurement in biological samples such as plasma or serum thus represents a suitable **surrogate marker** for novel Arginase inhibitors.

ASSAY SPECIFICATIONS

Format	96-well kit
Species Reactivity	Any species
Samples	Plasma, Serum, Cell culture
Sample volume	20 μ L
Sensitivity	2.1 μ M
Assay range	5.12 - 200 μ M
Assay time	Sample preparation: 3h ELISA overnight

STANDARD CURVE

Standard curve obtained with the hArg ELISA kit. In this competitive enzyme Immunoassay, optical density is inversely correlated with L-Arg levels within a linear range of 5.12 - 200 μ M.



METHOD VALIDATION

L-Arg was quantified in human plasma samples from 40 healthy donors either using IS-I-0400 ELISA kit or by liquid chromatography-mass spectrometry (LC/MS). Correlation study showed a $R^2 = 0.9713$, thereby confirming the accuracy of the immunoassay.

