

sHLA-G ELISA kit

The sHLA-G ELISA kit is a complete assay for the quantitative determination of soluble human HLA-G (leukocyte antigen-G) levels. This kit is to be used with EDTA plasma, amniotic fluid, and cell culture supernatant from human origin. The sHLA-G ELISA has strong cross-reactivity with bovine, dog, goat, monkey, rabbit and sheep HLA-G and weak cross-reactivity with cat HLA-G.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-309-KI01	96 wells
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

Application Notes For the quantitative determination of soluble human HLA-G (leukocyte antigen-G) in EDTA plasma, amniotic fluid, and cell culture supernatant.

Assay Time ~20 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Crossreactivity Cross-reacts strongly with bovine, dog, goat, monkey, rabbit and sheep HLA-G. Cross-reacts weakly with cat HLA-G.

Quantity For 96 wells (~80 tests)

Sensitivity 0.6U/ml (range 3.91 to 125U/ml)

Species Reactivity Human

Wavelength 450 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)