

AlphaLISA®

**Human IgG Fc Detection Kit**

**Product number:** AL3129 HV **Lot Number:** 3271365

---

**Material provided:****Kit Format:**

AL3129HV: 100 assay points

The number of assay points based on an assay volume of 100 µL in a 96-well assay plate using kit components at the recommended concentrations.

AL3129C: 500 assay points AL3129F: 5000 assay points

The number of assay points is based on an assay volume of 50 µL in 384-well assay plates using kit components at the recommended concentrations.

**Analyte Format:**

AL3129S: 200 µg (one vial contains an amount sufficient for performing 10 AlphaLISA standard curves)

**Manufacturing date:**

12/05/2023

**Document version:**

1

**Product Information****Application:**

This kit is designed for the quantitative determination of hIgG Fc in cell culture media and serum using a homogeneous AlphaLISA assay (no wash steps).

**Kit contents:**

The kit contains 5 components: AlphaLISA Acceptor beads coated with anti-hIgG Fc Antibody, Streptavidin-coated Donor beads, Biotinylated Anti-hIgG Fc Antibody, Lyophilized human IgG Fc and 10X AlphaLISA HiBlock Buffer.

**Sensitivity:**

Lower Detection Limit (LDL): 6 ng/mL

Lower Limit of Quantification (LLOQ): 35 ng/mL

EC<sub>50</sub>: 42 ng/mL**Dynamic Range:**

6 – 100 000 ng/mL

**Storage:**

Store kit in the dark at +4°C. The aliquoted analyte stored at -20 °C is stable up to 90 days. Avoid freezing analyte and freeze-thaw cycles.

**Stability:**

This kit is stable for at least 12 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

**Quality Control**

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals, EC<sub>50</sub> and LDL were measured on the EnVision Multilabel Plate Reader with Alpha option. We certify that these results meet our quality release criteria. Maximum counts may vary between bead lots and the instrument used, with no impact on LDL measurement.

EC <sub>50</sub> :	18.09 ng/mL
LDL:	6.19 ng/mL
LLOQ:	56.6 ng/mL
Min counts:	88 counts
Max counts:	157710 counts

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

AL3129-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000 [www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.