

AlphaLISA®

Human Immunoglobulin G subclass 1 (IgG₁ isotyping) Detection Kit**Product number:** AL3171 HV **Lot Number:** 3262931**Material provided:****Kit Format:**

AL3171HV: 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.

AL3171C: 500 assay points AL3171F: 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:

AL3171S: 10 µg (one vial contains an amount sufficient for performing 10 AlphaLISA standard curves)

Manufacturing date: February 14, 2024 **Document version:** 1**Product Information****Application:**

This kit is designed for the quantitative determination of human IgG₁ in cell culture/non-human serum, using a homogeneous AlphaLISA assay (no wash steps). The assay shows negligible cross-reactivity with other human IgG isotypes and monkey IgG.

Kit contents:

The kit contains 5 components: AlphaLISA Acceptor beads coated with anti-IgG₁ Antibody, Streptavidin-coated Donor beads, Biotinylated anti-IgG₁ antibody, Lyophilized Human IgG₁, and 10X AlphaLISA HiBlock Buffer.

Sensitivity:

Lower Detection Limit (LDL): 3.82 ng/mL

Lower Limit of Quantification (LLOQ): 4.29 ng/mL

EC₅₀: 3.68 µg/mL**Dynamic Range:**

4.29 – 1 000 000 ng/mL

Storage:

Store kit in the dark at +4°C. Store reconstituted analyte at -20°C.

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. Note: Once reconstituted, the human IgG₁ analyte is stable for at least 1 months when stored at -20°C.

Quality Control

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals, EC₅₀ 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 ₅₀ :	3.218 µg/mL
LDL:	4.006 ng/mL
LLOQ:	11.987 ng/mL
Min counts:	105 counts
Max counts:	129900 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.

AL3171-R Rev01

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a modern, clean appearance.

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

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.