

AlphaLISA®

Human Corticosterone Detection Kit

Product number: AL3020 HV **Lot Number:** 3223067

Material provided:

Kit Format: AL3020HV: 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.

AL3020C: 500 assay points AL3020F: 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: AL3020S: 100 µL at 10 mg/mL (one vial contains an amount sufficient for performing 8 AlphaLISA standard curves)

Manufacturing date: December 8, 2023 **Document version:** 1

Product Information

Application: This kit is designed for the quantitative determination of Corticosterone in serum and plasma of mammalian species and chickens using a homogeneous AlphaLISA assay (no wash steps). The assay shows minimum cross reactivity with other steroids.

Kit contents: The kit contains 5 components: AlphaLISA Acceptor beads coated with Anti-CORT antibody, Streptavidin-coated Donor beads, Biotinylated CORT-BSA Tracer, CORT Analyte, and 5X NaCl Buffer.

Sensitivity: Lower Detection Limit (LDL): 9 ng/mL
IC₅₀: 37 ng/mL

Dynamic Range: 0.02 – 8000 ng/mL

Storage: Store kit in the dark at +4°C except the analyte that should be stored at room temperature. It has been demonstrated that CORT DMSO solution is stable for at least 6 months at room temperature.

Stability: This kit is stable for at least 6 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, 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.

IC ₅₀ :	25.130 ng/mL
LDL:	10.121 ng/mL
Min counts:	170 counts
Max counts:	25920 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.

AL3020-R Rev01

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a slightly irregular, hand-drawn 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.