

AlphaLISA®

Di-Methyl-Histone H3 Lysine 36 (H3K36me2) Cellular Detection Kit**Product number:** AL723 F **Lot Number:** 3281595**Material provided:****Kit Format:**

AL723HV: 100 assay points

The number of assay points based on an assay volume of 50 µL in a 96-well 1/2 area assay plate using kit components at the recommended concentrations.

AL723C: 500 assay points AL723F: 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.

Manufacturing date: March 29, 2024 **Document version:** 1**Product Information****Kit contents:**

The kit contains 6 components: AlphaLISA Acceptor beads coated with an anti-epigenetic mark antibody, Streptavidin-coated Donor beads, Biotinylated anti-Histone H3 (C-terminus) Antibody, and Cell-Histone™ Lysis (1X), Extraction (1X) and Detection (10X) buffers.

Storage:

Store kit in the dark at 4 °C.

Stability:

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

Application:

This kit is designed for the detection of di-methylated Histone H3 Lysine 36 (H3K36me2) in cell lysates using a homogeneous AlphaLISA assay (no wash steps).

Quality Control

Lot-to-lot consistency of Donor and Acceptor beads is confirmed by a Quality Control AlphaLISA titration assay read on an EnVision® instrument. Maximum signal and EC₅₀ value are determined using a biotin-H3K36me2 peptide. Minimum signal is measured in the absence of peptide. Maximum counts may vary between bead lots. Maximum counts obtained in the QC assay are usually higher than those obtained in a cellular detection assay, which are dependent on epigenetic mark abundance and assay conditions (e.g. cell line, culture medium, incubation time, modulator concentration, etc.).

Maximum signal: 1,115,572 counts
Minimum signal: 348 counts
EC₅₀: 8.18 xM

QC release specifications of the biotinylated antibody are based on spectrophotometric analysis of the labeled antibody.

Labeling Ratio: 4.56 biotin/Ab

We certify that these results meet our quality release criteria.

Please visit our website for additional information on AlphaLISA technology at www.revivity.com

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.

AL723-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.revivity.com

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