

AlphaLISA®

Epigenetics Buffer 1 (5X)

Product number: AL008F

Lot Number: 3345752

Product Format:

Kit Components	10 mL Kit	100 mL Kit
AlphaLISA 5X Epigenetics Buffer 1	AL008C1 (10 mL)	AL008F1 (100 mL)
AlphaLISA 30X Epigenetics Buffer Supplement	AL008C2 (1.7 mL)	AL008F2 (17 mL)

Manufacturing date: October 15, 2024

Document version: 1

Product Information

Description: This product is a two-component buffer kit specifically formulated for the detection of histone H3-derived peptides in AlphaLISA Epigenetics assays.

Application: Once prepared, the 1X AlphaLISA Epigenetics Buffer 1 is used for diluting AlphaLISA Acceptor and Alpha Donor beads. It prevents the non-specific binding of highly-charged histone peptides to the Acceptor and Alpha Donor beads in AlphaLISA Epigenetics assays.

Storage: Store kit components at 4°C.

Stability: The undiluted stock solutions are stable for at least 12 months from the manufacturing date when stored in their original packaging at the recommended storage conditions. Upon preparation, the 1X AlphaLISA Epigenetics Buffer 1 should be used within 16 hours.

Preparation: To 7.67 volumes of ultrapure water (18 MΩ•cm), add 2 volumes of 5X Epigenetics Buffer 1 and 0.33 volume of 30X Buffer Supplement. The 1X AlphaLISA Epigenetics Buffer 1 may appear slightly cloudy or precipitation may occur. However, this will not affect assay performance. DO NOT heat above 30°C as this could significantly affect performance.

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

AL008-R Rev 01