

AlphaLISA®

Bovine Immunoassay Buffer (10X)

Product number: AL010C

Lot Number: 3382052

Product Format: AL010C: 10 mL

AL010F: 100 mL

This product is designed to be used in combination with Bovine IgG, IgG1, IgG2, IgA, IgM, IFN γ , and TNF α AlphaLISA kits.
Manufacturing date: February 10, 2025

Document version: 1

Product Information

Description: AlphaLISA Bovine Immunoassay Buffer (10X)

Presentation: Liquid, 10X

Formulation (10X): 10X PBS, pH 7.3 – 7.4, 1.0% Tween-20, and 0.5% Kathon as a preservative

Storage: Store 10X AlphaLISA Bovine Immunoassay Buffer at 4°C

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

Reconstitution: Add 1 volume of 10X AlphaLISA Bovine Immunoassay Buffer to 9 volumes of Milli-Q® H2O. 1X Buffer is stable for 30 days at 4°C. Long-term storage of 1X Buffer is not recommended.

Once diluted, 1X AlphaLISA Bovine Immunoassay Buffer contains 1X PBS, pH 7.4, 0.1% Tween-20 and 0.05% Kathon.

 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.

AL010-R Rev 01