

AlphaLISA®

## ImmunoAssay Buffer (10X)

**Product number:** AL000C **Lot Number:** 3198914

**Product Format:** AL000C: 10 mL  
AL000F: 100 mL

**Manufacturing date:** 9/18/2023 **Document version:** 1

### Product Information

**Description:** AlphaLISA ImmunoAssay Buffer (10X)

**Presentation:** Liquid, 10X  
*10X buffer might be slightly yellow. However, this does not affect the assay results.*

**Formulation (10X):** 250 mM HEPES, pH 7.4, 1% Casein, 10 mg/mL Dextran-500, 5% Triton X-100 and 0.5% Kathon

**Storage:** Store 10X AlphaLISA ImmunoAssay Buffer at 4°C

**Stability:** This product is stable for at least 24 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

**Reconstitution:** Add 1 volume of 10X AlphaLISA ImmunoAssay Buffer to 9 volumes of dd H<sub>2</sub>O or Milli-Q® H<sub>2</sub>O. 1X Buffer is stable for 3 days at 4°C. Long-term storage of 1X Buffer is not recommended.

Once diluted, 1X AlphaLISA ImmunoAssay Buffer contains 25 mM HEPES, pH 7.4, 0.1% Casein, 1 mg/mL Dextran-500, 0.5% Triton X-100 and 0.05% Kathon.

Please visit our website for additional information on AlphaLISA technology at [www.revvity.com](http://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.

AL000-R Rev 01