

AlphaLISA®

HiBlock Buffer (10X)

Product number: AL004C Lot Number: 3321716

Product Format: AL004C: 10 mL
AL004F: 100 mL

Manufacturing date: August 1, 2024 Document version: 1

Product Information

- Application:** This product is designed to be used in combination with AlphaLISA kits. HiBlock buffer was especially developed buffer to block non-specific protein binding sites.
- Formulation (10X):** 250 mM HEPES, pH 7.4, 1% Casein, 10 mg/mL Dextran-500, 5% Triton X-100, 5% Blocking reagent, 5% BSA and 0.5% Kathon.
- Storage:** Store 10X AlphaLISA HiBlock 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. Once diluted, 1X AlphaLISA HiBlock Buffer is stable for 3 days at +4°C, long term storage of 1X buffer is not recommended.
- Reconstitution:** Add 1 volume of 10X AlphaLISA HiBlock Buffer to 9 volumes of dd H₂O or Milli-Q® H₂O.
Once diluted, 1X AlphaLISA HiBlock Buffer contains: 25 mM HEPES, pH 7.4, 0.1% Casein, 1 mg/mL Dextran-500, 0.5% Triton X-100, 0.5% Blocking reagent, 0.5 % BSA and 0.05% Kathon.
- Appearance:** 10X buffer is slightly yellow-brownish (Figure 1).



Fig.1. Typical appearance of 10X HiBlock Buffer.

Notes

- Occasionally, turbidity can occur in HiBlock buffer. This cloudy appearance is caused by flocculation or coagulation of components present in the buffer (Figure 2), especially at cold temperatures, which could occur during shipping.

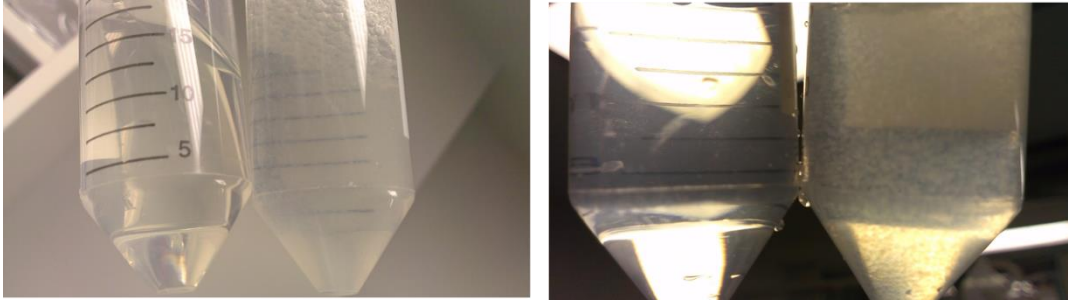


Fig.2.

HiBlock buffer when fully dissolved (left tube on each image) and containing gelatin in suspension (right tube on each image).

- Should flocculation or precipitation of suspensions of buffer components occur, it is recommended to dilute to a 1x solution and centrifuge at 1000 rpm for 5 min to pellet the precipitates.
- The appearance of the buffer does not affect its efficacy (Figure 3).

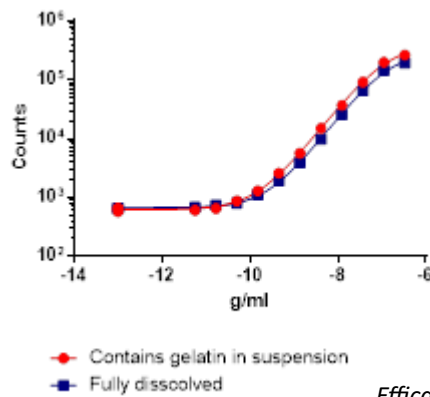


Fig. 3.

Shown are results generated with the Human IgG4 kit.

Efficacy of clear versus turbid HiBlock buffer.

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.

AL004-R Rev01

revvity