

HiBlock Buffer

Caution: For Laboratory Use. A research chemical for research purposes only.

Product information

Application: This product is designed to be used in combination with AlphaLISA kits. HiBlock Buffer was especially developed to block non-specific protein binding sites.

Formulation: 10X AlphaLISA HiBlock Buffer contains: 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% Proclin-300.

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% Proclin-300.

Storage: Store 10X HiBlock Buffer at +4 °C. 1X HiBlock Buffer is stable for 3 days at +4 °C.

Appearance: 10X HiBlock Buffer is slightly yellow-brownish (Figure 1) in appearance.

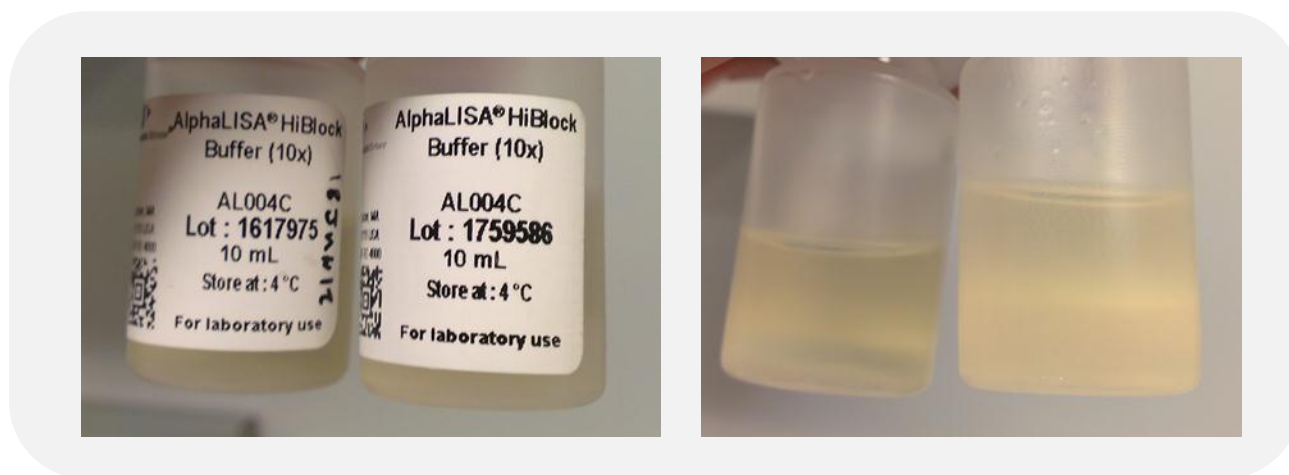


Figure 1. Typical appearance of 10X HiBlock Buffer.

Notes

- Flocculation or precipitation can occur during shipping (Figure 2), especially at cold temperatures.

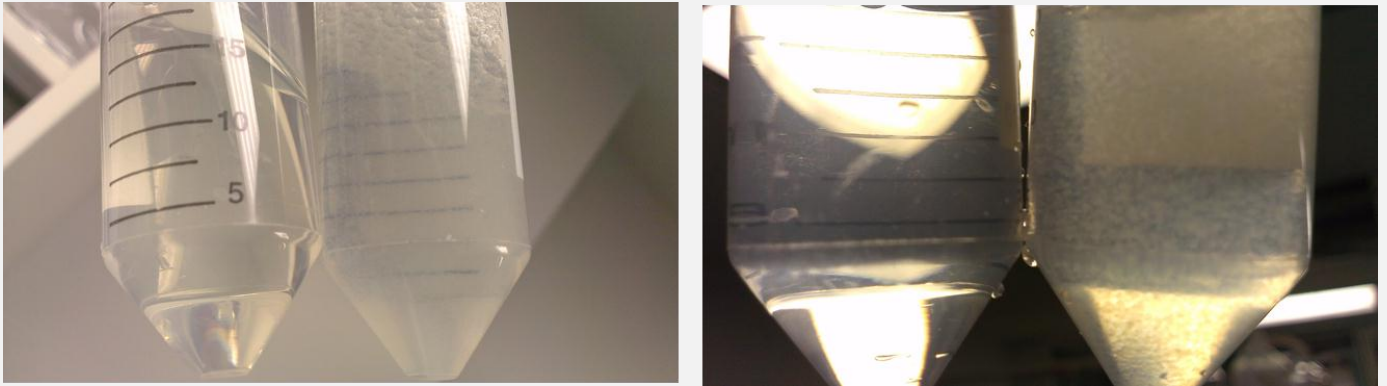


Figure 2. HiBlock Buffer when fully dissolved (left tube on each image) and containing caseins in suspension (right tube on each image).

- The appearance of the buffer does not affect its efficacy (Figure 3).

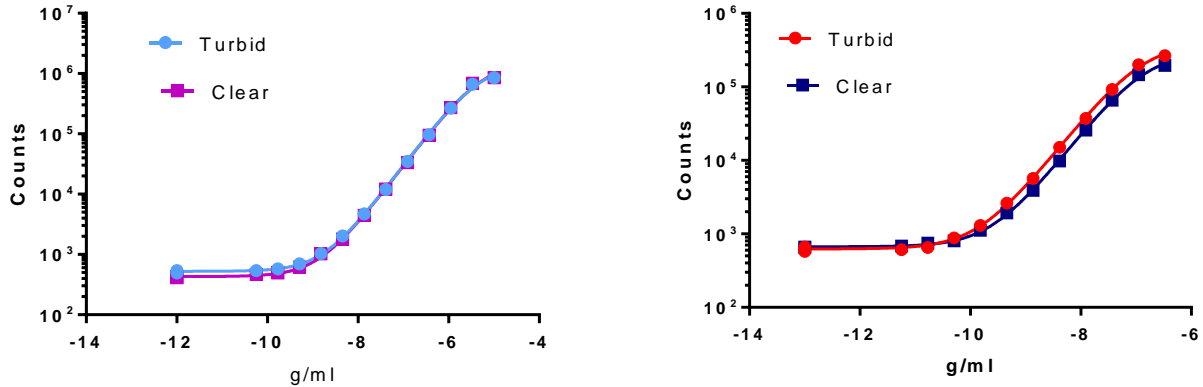


Figure 3. Efficacy of clear versus turbid HiBlock Buffer using the human Transferrin AlphaLISA kit (left) and human IgG4 kit (right).

- However, 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.
- Warming to 37°C will help solubilize the suspension without completely removing precipitates.

A comparative study was performed using the Transferrin AlphaLISA kit in the presence of turbid HiBlock Buffer used as is, or after heating (Warm) or centrifugation (SpinDown). As shown in Figure 4, all 3 conditions yielded identical results although the appearance of the centrifuged buffer was best.

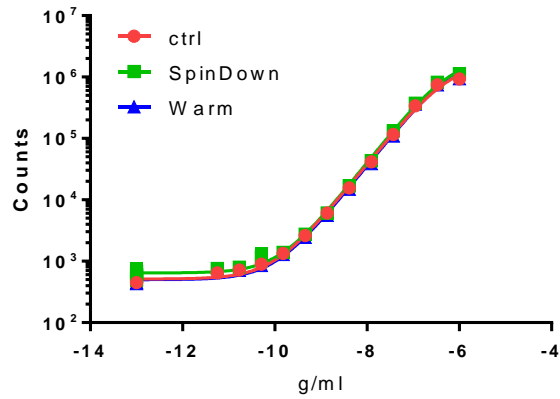


Figure 4. Effect of different treatments on the efficacy of HiBlock Buffer in the Transferrin AlphaLISA assay.

Please visit our website for more information:

<http://www.perkinelmer.com/in/Catalog/Family/ID/AlphaLISA%20HiBlock%20Buffer>

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. Corning and Epic are registered trademarks of Corning Incorporated. All other trademarks are the property of their respective owners.