

LANCE® *Ultra***HiBlock Buffer (5X)**

**Product number:** TRF1011C **Lot Number:** 3262933

**Product Format:** TRF1011C: 10 mL  
TRF1011F: 100 mL

**Manufacturing date:** February 8, 2024 **Document version:** 1

**Product Information**

- Application:** This product is designed to be used in combination with LANCE *Ultra* kits. The *Ultra* HiBlock buffer was especially developed buffer to block non-specific protein binding sites.
- Formulation (5X):** 5X *Ultra* HiBlock Buffer contains: 125 mM HEPES, pH 7.4, 0.5% Casein, 5 mg/mL Dextran-500 K, 2.5% Triton X-100, 2.5% Blocking reagent, 2.5% BSA and 0.25% Kathon.
- Storage:** Store 5X Buffer at 4°C. 1X *Ultra* HiBlock Buffer is stable for 30 days at 4°C. Long-term storage of 1X *Ultra* HiBlock Buffer is not recommended.
- 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, *Ultra* Buffer is stable for 3 days at +4°C, long term storage of 1X buffer is not recommended.
- Reconstitution:** Add 2 volumes of 5X *Ultra* HiBlock Buffer to 8 volumes of Milli-Q® H<sub>2</sub>O.  
Once diluted, 1X *Ultra* HiBlock Buffer contains: 25 mM HEPES, pH 7.4, 0.1% Casein, 1 mg/mL Dextran-500K, 0.5% Triton X-100, 0.5% Blocking reagent, 0.5 % BSA and 0.05% Kathon.
- Appearance:** 5X *Ultra* HiBlock Buffer is slightly tan (Figure 1).



Fig.1. Typical appearance of 5X *Ultra* HiBlock Buffer.

Please visit our website for additional information on LANCE Ultra technology at [www.revivity.com](http://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.

TRF1011-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000 [www.revivity.com](http://www.revivity.com)

For a complete listing of our global offices, visit [www.revivity.com](http://www.revivity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.