

HTRF® Research Reagents

HTRF Epig. Detect. Buf. 1 - 24 mL

Product number: 62SAHRDD

Lot number: 15RA

Product characteristics:

- Conditioning buffer -
- Appearance Frozen form
- Storage conditions <-16 °C
- Expiration date 2024.10.09
- Volume per vial 24.00 mL

Weight indications:

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (Buf) | 24.00 mL | N/A |

Please note

The expiration date of batch 15RA has been extended to 2024-10-09 after validation by QC.

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.