

HTRF® Research Reagents

HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL

Product number: 64KL1FDF

Lot number: 34RA

Product characteristics:

• Conditioning buffer	Stock solution 4X
• Appearance	Liquid form
• Storage conditions	2-8 °C
• Expiration date	2027.12.15
• Volume per vial	130.00 mL

Weight indications:

	Quantity per vial	MW (Da)
Active Moiety (Buf)	130.00 mL	N/A

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.