

HTRF® Research Reagents

HTRF Europium Cryptate Labeling kit

Product Number : 62EUSPEA Lot Number : 27A

Expiration date : 2026-10-20 Document version : 1

Reagents :	Lot n°
HTRF Eu cryptate - 620 nm calibrator	19RA
HTRF Purification Column 1	45A
HTRF Elution buf. 1 - 20 mL	16RA

HTRF Europium Cryptate Labeling kit - Lot n° : 27A

 Accepted

 Rejected

Quality control date : 2024-09-02	Supervisor's initials : MP
-----------------------------------	----------------------------

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.