

HTRF® Research Reagents

HTRF MAb Anti p-AKT-Eu - 500 pts

Product number: 61P15KAE

Lot number: 08A

Product characteristics:

• Packaging Unit	500 tests/vial
• Conditioning buffer	Hepes 0.05M pH7 + 0.1% BSA
• Appearance	Frozen form
• Storage conditions	-20 °C
• Expiration date	2027.07.24
• Volume per vial	250.00 µL
• Active Moiety Concentration (Ab)	6.00 µg/mL

Weight indications:

	Quantity per vial	MW (Da)
Active Moiety (Ab)	1.50 µg	150 000

Recommended quantity per well*:

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	3.00 ng
Number of tests/vial	500

*Actual amount per well should be optimized according to assay conditions

Please note

The expiration date of batch 08A has been extended to 2027-07-24 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.