

HTRF® Research Reagents

HTRF PAb Anti Human IgG-Tb - 5K pts
Product number: 61HFCTAA

Lot number: 05A

Product characteristics:

- | | |
|--|-------------------------------|
| • Packaging Unit | 5 000 tests/vial |
| • Conditioning buffer | Phosphate 0.1M pH7 + 0.1% BSA |
| • Appearance | Frozen form |
| • Storage conditions | <-16 °C |
| • Expiration date | 2033.05.08 |
| • Volume per vial | 0.25 mL |
| • Active Moiety Concentration (PAb) | 40.00 µg/mL |

Weight indications:

	Quantity per vial	MW (Da)
Active Moiety (PAb)	10.00 µg	150 000

Recommended quantity per well*:

	in 384 well low volume plate (20µL)
Active Moiety (PAb)	2.00 ng
Number of tests/vial	5 000

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

Please note

The use of KF is not compulsory in HTRF assays using Lumi4-Tb conjugate.

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.