

**phosph p53-K**  
Product reference :

anti phospho P53 cryptate conjugate  
**61P08KAZ**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 05
- Packaging Unit\_\_\_\_\_ 10 000 tests/vial
- Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.1% BSA + 0.1% tween
- Appearance\_\_\_\_\_ Frozen form
- Storage conditions\_\_\_\_\_ <-16 °C
- Expiration date\_\_\_\_\_ 2024.03.02
- Volume per vial\_\_\_\_\_ 250 µL
- Active Moiety Concentration (Ab)\_\_\_\_\_ 100 µg/mL

**Weight indications :**

|                    | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (Ab) | 25.00 µg          | 150 000 |

**Recommended quantity per well\* :**

|                      | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (Ab)   | 2.50 ng                             |
| Number of tests/vial | 10 000                              |

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

**PLEASE NOTE**

KF can play an essential role in the lanthanide cryptate protection by preventing the action of possible quenchers contained in the assay. KF is mandatory for HTRF assays using Europium cryptate. KF is generally used at a final concentration of 100 to 400 mM, and is added to the conjugate working solutions or dispensed in a separate step, just before the readout.

**Further information as well as a specific application note regarding conjugate-related calculations are available on our HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)**