

**UBIQ-K**

Ubiquitin-Eu cryptate - 5,000 tests

Product reference :

**61UBIKLA****Product characteristics :**

- Lot Number	11A
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.5% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2022.02.08
- Reconstitute each vial with	0.25 mL distilled H2O

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ubq)	3.75 µg	8 500

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Ubq)	751.00 pg
Number of tests/vial	5 000

\*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.

The expiration date of batch 11A has been extended to 2022-02-08 after validation by QC

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