

Batch information

Mab HA-K

HTRF MAb Anti HA-Eu - 20K pts

Product reference :

610HAKLB

Product characteristics :

- Lot Number	14A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.5% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2023.11.05
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (Mab)	16.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	16.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	0.80 ng
Number of tests/vial	20 000

*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 09Z has been extended to 2021-10-14 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

