

## Batch information

### Anti-GST-K

MAB Anti GST-Eu cryptate - 5,000 tests

Product reference :

**61GSTKLA**

### Product characteristics :

- Lot Number	104A
- Conditioning buffer	Phosphate 0.1M pH7 + 0.5% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.03.03
- Reconstitute each vial with	250 µL distilled H2O
- Active Moiety Concentration (Mab)	50.00 µg/mL

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	12.50 µg	150 000

### 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 Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)**

