

Anti-GST-K

MAb Anti GST-Eu cryptate - 5,000 tests

Product reference :

61GSTKLA**Product characteristics :**

- Lot Number_____	101B
- Conditioning buffer_____	Phosphate 0.1M pH7 + 0.5% BSA + stabilizers
- Appearance_____	Lyophilized form
- Storage conditions_____	2-8 °C
- Expiration date_____	2023.01.29
- Reconstitute each vial with_____	250 µL distilled H2O
- Active Moiety Concentration (Mab)_____	50 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. The expiration date of batch 101A has been extended to 2022-05-10 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