

Anti-GST-K

MAb Anti GST-Eu cryptate - 1 mg

Product reference :

61GSTKAH

Product characteristics :

- Lot Number_____	100A
- Packaging Unit_____	400 000 tests/vial
- Conditioning buffer_____	Phosphate 0.1M pH7 + 0.1% BSA
- Appearance_____	Frozen form
- Storage conditions_____	<-16 °C
- Expiration date_____	2025.08.16
- Reconstitute each vial with_____	4 mL distilled H2O
- Volume per vial_____	4 mL
- Active Moiety Concentration (Mab)_____	250 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	1.00 mg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	2.50 ng
Number of tests/vial	400 000

*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