

61KAPKAB

MAb Anti Human Kappa-Eu cryptate - 20,000 tests - CUSTOM LABELING

Product reference :

61KAPKAB**Product characteristics :**

- Lot Number	006
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	-80 °C
- Expiration date	2028.09.28
- Volume per vial	200 µL
- Active Moiety Concentration (Ab)	100 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	20 000.00 ng	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	1.00 ng
Number of tests/vial	20 000

*Actual amount per well should be optimized according to assay conditions

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com