

61KAPDAA

MAb Anti Human Kappa-d2 - 5,000 tests - CUSTOM LABELING

Product reference :

61KAPDAA**Product characteristics :**

- Lot Number	006
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	-80 °C
- Expiration date	2029.01.13
- Volume per vial	50 µL
- Active Moiety Concentration (Ab)	500 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	25 000.00 ng	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	5.00 ng
Number of tests/vial	5 000

*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