

Batch information

Anti-Mouse - d2

HTRF PAb Anti Mouse IgG-d2 - 1K pts

Product reference :

61PAMDAF

Product characteristics :

- Lot Number	03B
- Packaging Unit	1 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2030.10.11
- Volume per vial	0.08 mL
- Active Moiety Concentration (PAb)	250.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (PAb)	20.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (PAb)	20.00 ng
Number of tests/vial	1 000

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

PLEASE NOTE

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

