

Batch information

Anti-Mouse-Tb

HTRF PAb Anti Mouse IgG-Tb - 5K pts

Product reference :

61PAMTAA

Product characteristics :

- Lot Number	03B
- Packaging Unit	5 000 tests/vial
- Conditioning buffer	Hepes 0.05 M pH7 + 0.1 % BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2024.11.09
- Volume per vial	0.25 mL
- Active Moiety Concentration (Ab)	25.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	6.25 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	1.25 ng
Number of tests/vial	5 000

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

PLEASE NOTE

The use of KF is not compulsory in HTRF assays using Lumi4-Tb conjugate.

The expiration date of batch 03B has been extended to 2024-11-09 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

