

PAb Anti human IgG-Tb
Product reference :

PAb Anti Human IgG-Tb cryptate - 1,000 tests
61HFCTAF

Product characteristics :

- Lot Number _____ 02A
- Packaging Unit _____ 1 000 tests/vial
- Conditioning buffer _____ Phosphate 0.1M pH7 + 0.1% BSA
- Appearance _____ Frozen form
- Storage conditions _____ <-16 °C
- Expiration date _____ 2028.02.05
- Volume per vial _____ 50 µL
- Active Moiety Concentration (PAb) _____ 40 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (PAb)	2.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (PAb)	2.00 ng
Number of tests/vial	1 000

*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.

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