

**PAb Anti human IgG-Tb**

Product reference :

PAb Anti Human IgG-Tb cryptate - 1 mg

**61HFCTAH**

**Product characteristics :**

- Lot Number	01A
- Packaging Unit	500 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% BSA
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2025.07.23
- Volume per vial	4 mL
- Active Moiety Concentration (PAb)	250 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (PAb)	1.00 mg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (PAb)	2.00 ng
Number of tests/vial	500 000

\*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 HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)**