

MAB Anti 6His Tb cryptate Gold MAb Anti-6HIS Tb cryptate Gold - 1 mg - CUSTOM LABELING
 Product reference : **61HI2TAH**

Product characteristics :

- Lot Number_____ 03A
- Conditioning buffer_____ P04 100 mM pH7 + BSA 0.1% + Tween 0.1%
- Appearance_____ Frozen form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2026.04.13
- Volume per vial_____ 3.402 mL
- Active Moiety Concentration (HIS)_____ 294 µg/mL

Weight indications :

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

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

