

Lumi 4 Tb-NHS
Product reference :

Terbium cryptate NHS - 0.1 mg
65TBSABA

Product characteristics :

- Lot Number_____ 26A
- Conditioning buffer_____ DMSO
- Appearance_____ Frozen form
- Storage conditions_____ <-60 °C
- Expiration date_____ 2023.04.18
- Volume per vial_____ 10.67 µL
- Active Moiety Concentration (K)_____ 6.28 mM

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (K)	0.10 mg	1 529

PLEASE NOTE

0.1 mg corresponds to 67 nmoles of Lumi 4 Tb-NHS
Rz (337/280) = 7.7 Molar extinction coefficient (337nm) = 26 000 L.Mole⁻¹
Caution : this material is for internal use only !

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