

KTBP5COOHMP

Europium cryptate TBP 5COOH MP - 0.1 mg

Product reference :

65T5MABA**Product characteristics :**

- Lot Number _____ 07A
- Conditioning buffer _____ Desiccated
- Appearance _____ Desiccated form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2022.09.12

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (K)	0.10 mg	1 523

PLEASE NOTE

Rz (320.5/280) = 2.82

Molar extinction coefficient (317 nm) = 31 000 L.Mole⁻¹

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