

S0-500 biotin
Product reference :

TK Substrate-biotin - 500 µg
61TK0BLC

Product characteristics :

- Lot Number	16A
- Conditioning buffer	Hepes buf. 20mM PH7.4, BSA 1%+ stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2022.06.19

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (sub)	500.00 µg	1 740

PLEASE NOTE

Reconstitute with 574µL H₂O to obtain a 500µM stock solution -Refer to kinEASE TK bulk HTRF package insert (ref .doc. 62TK0PEC) for more information

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