

S0-500 biotin
Product reference :

TK Substrate-biotin - 500 µg
61TK0BLC

Product characteristics :

- Lot Number_____ 21A
- Conditioning buffer_____ Hepes buf. 20mM PH7.4,
BSA 1%+ stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2023.10.08

Weight indications :

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

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