

S1-Biotin

STK-Substrate 1-biotin - 50 µg

Product reference :

61ST1BLE

Product characteristics :

- Lot Number _____ 07A
- Conditioning buffer _____ Hépes buffer 20 mM pH 7.4 +
BSA 1% + stabilizers
- Appearance _____ Lyophilized form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2022.09.11

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Sub)	50.00 µg	1 687

PLEASE NOTE

Reconstitute with 593µL H₂O to obtain a 50µM stock solution -Refer to kinEASE HTRF package insert (ref .doc. 62STxPEB) 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

