

Batch information

STD HTRF BDNF kit - Std
Product reference : **62BDNFCDA**

Product characteristics :

| | |
|------------------------------------|--|
| - Lot Number | 01A |
| - Conditioning buffer | Phosphate 50mM pH7.0 +BSA 1% + stabilizers |
| - Appearance | Lyophilized form |
| - Storage conditions | 2-8 °C |
| - Expiration date | 2024.09.01 |
| - Active Moiety Concentration (Ab) | 15.00 ng/mL |

Weight indications :

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (Ab) | 4.50 ng | N/A |

PLEASE NOTE

Reconstitute with 300 µl H₂O

The expiration date of batch 01A has been extended to 2024-09-01 after validation by QC.

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

