

**STD**

BDNF kit - Standard

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\_\_\_\_\_ 2022.09.01  
- Active Moiety Concentration (Ab)\_\_\_\_\_ 15 ng/mL

**Weight indications :**

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

**PLEASE NOTE**

Reconstitute with 300 µl H<sub>2</sub>O

***Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)***

