

**Mouse STING WT bind kit - Std**      Mouse STING WT binding kit - Standard  
Product reference :                      **64BDSTGMCA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 01A  
- Conditioning buffer \_\_\_\_\_ Hepes buffer pH7 - BSA -  
preservatives  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ <-60 °C  
- Expiration date \_\_\_\_\_ 2022.08.23  
- Volume per vial \_\_\_\_\_ 50 µL  
- Active Moiety Concentration (Pro) \_\_\_\_\_ 2 µM

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Pro)	0.10 nmoles	N/A

***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)***