

**STD**

Human STING WT binding kit - Standard  
**64BDSTGCDA**

Product reference :

**Product characteristics :**

- Lot Number \_\_\_\_\_ 03A  
- Conditioning buffer \_\_\_\_\_ Tampon hépés + stabilizer  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ -60 °C  
- Expiration date \_\_\_\_\_ 2023.01.02  
- Volume per vial \_\_\_\_\_ 50 µL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	10.00 nmoles	N/A

**PLEASE NOTE**

*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)*

