

**mTOR phospho-S2448 Kit control lysate** / **mTOR phospho-S2448 kit control lysate**  
Product reference : **64TORTDA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 05A  
- Conditioning buffer \_\_\_\_\_ Tampon de lyse compl ment  (1X)  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ -60  C  
- Expiration date \_\_\_\_\_ 2022.02.24

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiey (Ab)	150.00 �l	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)*