

**Control lysate**

FoxO1 phospho-S256 kit control lysate

Product reference :

**64FOXTDA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 02A  
- Conditioning buffer \_\_\_\_\_ Lysis Buffer  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ -60 °C  
- Expiration date \_\_\_\_\_ 2022.12.03  
- Volume per vial \_\_\_\_\_ 150 µL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	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)*