

**Control lysate**

AMPK total kit control lysate

Product reference :

**63ADK060TDA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 07A  
- Conditioning buffer \_\_\_\_\_ lysis buffer  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ -60 °C  
- Expiration date \_\_\_\_\_ 2022.09.04  
- 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)*