

**Phospho ACC control lysate**

ACC phospho-S79 kit - control lysate

Product reference :

**64ACCTDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 12A  
- Conditioning buffer\_\_\_\_\_ lysis buffer  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ -60 °C  
- Expiration date\_\_\_\_\_ 2023.09.06  
- Volume per vial\_\_\_\_\_ 150 µL  
- Active Moiety Concentration (PHO)\_\_\_\_\_ 1 X

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (PHO)	150.00 µl	N/A

**PLEASE NOTE**

The expiration date of batch 12A has been extended to 2023-09-06 after validation by QC

***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)***