

**Phospho ACC control lysate**

ACC phospho-S79 kit - control lysate

Product reference :

**64ACCTDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 11A  
- Conditioning buffer\_\_\_\_\_ lysis buffer  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ -60 °C  
- Expiration date\_\_\_\_\_ 2023.04.03  
- 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 11A has been extended to 2023-04-03 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)***