

**ERK phospho-T202/Y204 kit control lysate** ERK phospho-T202/Y204 kit - control lysate  
Product reference : **62ERKTDA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 13A  
- Conditioning buffer \_\_\_\_\_ Phosphate 0.1M pH7 + % BSA +  
% Tween  
- Appearance \_\_\_\_\_ Lyophilized form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2023.09.10

**Weight indications :**

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

**PLEASE NOTE**

Reconstitute with 0.500 mL H<sub>2</sub>O.

***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)***