

## Batch information

ERK phospho-T202/Y204 kit control | ERK phospho-T202/Y204 kit - control lysate

Product reference : **62ERKTDA**

### Product characteristics :

- Lot Number	14A
- Conditioning buffer	Phosphate 0.1M pH7 + % BSA + % Tween
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2024.01.28

### Weight indications :

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

### PLEASE NOTE

Reconstitute with 0.400 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)*

