

## Batch information

### control lysate

HTRF ATF2 p-T71 kit - Ctrl lysate

Product reference :

**63ADK015TDA**

### Product characteristics :

- Lot Number	06A
- Conditioning buffer	Phosphate 0.1M pH7 + % BSA + % Tween
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2022.10.12

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	150 000

*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)*

