

## Batch information

**TYK2 total control lysate**

HTRF TYK2 Total Kit - Ctrl Lysate

Product reference :

**64TYK2TTDA**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	Lysis buffer 4
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.12.15

### Weight indications :

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

**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)**

