

Batch information

TYK2 p-Y1054/1055 control lysate HTRF TYK2 p-Y1054/1055 Kit - Ctrl Lysate

Product reference : **64TYK2Y10TDA**

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

