

Batch information

Control lysate

HTRF JAK2 p-Y1007/08 Kit - Ctrl lysate

Product reference :

64JAK2Y10TDA

Product characteristics :

- Lot Number	01A
- Conditioning buffer	Lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.09.22
- Volume per vial	150.00 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (CL)	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

