

Batch information

Control lysate HTRF JAK2 Total Kit - Ctrl lysate
Product reference : **64JAK2TTDA**

Product characteristics :

| | |
|-----------------------|--------------|
| - Lot Number | 01A |
| - Conditioning buffer | Lysis Buffer |
| - Appearance | Frozen form |
| - Storage conditions | -60 °C |
| - Expiration date | 2024.09.13 |
| - 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

