

Batch information

Control lysate

HTRF GAPDH House. kit - Ctrl lysate

Product reference :

64GAPDHTDA

Product characteristics :

| | |
|------------------------------------|--------------|
| - Lot Number | 03A |
| - Conditioning buffer | Lysis Buffer |
| - Appearance | Frozen form |
| - Storage conditions | -60 °C |
| - Expiration date | 2024.11.29 |
| - Volume per vial | 150.00 µL |
| - Active Moiety Concentration (CL) | 1.00 X |

Weight indications :

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (CL) | 150.00 µl | N/A |

PLEASE NOTE

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

