

**Control lysate**

STAT3 total kit control lysate

Product reference :

**64NT3TDA**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 07A  
- Conditioning buffer \_\_\_\_\_ Lysis Buffer  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ -60 °C  
- Expiration date \_\_\_\_\_ 2023.02.13  
- Volume per vial \_\_\_\_\_ 150 µL

**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)*

